

revvity

LabChip User Guide

ProteinEXact™ HR Assay User Guide

For LabChip® GXII Touch

©Copyright 2022-2024 Revvity Inc. All rights reserved. PN CLS150849 Rev G
Publication Date: June 10, 2024

CE



Table of Contents

Specifications	3
Assay Specifications	3
Sample Conditions	4
Storage Conditions	4
Reagent Kit Contents	5
LabChips	5
Safety and Usage	6
Safety Warnings and Precautions	6
Usage	6
Preparation Procedures	7
Additional Items Required	7
Preparing the Gel-Dye Solution	8
For High Throughput Chip Preparation (up to 96 samples)	8
For Low Throughput Chip Preparation (48 samples)	9
Preparing the Ladder and Wash Buffer Tube	10
Preparing the Chip	11
Inserting a Chip into the LabChip GXII Touch	13
Priming and Warming	14
Preparing the Protein Samples	16
Running the Assay	17
Cleaning and Storing the Chip	20
Chip Cartridge Cleaning	21
Results	22
ProteinEXact HR Ladder Result	22
Troubleshooting	23
LabChip Kit Essential Practices	30
General	30
Reverse Pipetting Technique	31
Reagents	32
Chips	33
General Guidelines	33
Repriming Chips	33
Washing Chips	34
Removing Sipper Clogs	36
Samples	36
Compatible Buffers, Salts and Additives	37
Chip Well Aspiration Using a Vacuum	38
Customer Technical Support	39
Licenses and Rights of Use	39

Specifications

Assay Specifications

Table 1. Assay Specifications

Sizing Range	6.5 - 250 kDa
Linear Concentration Range (R^2)	0.99 (10 - 1000 ng/ μ L) and 0.98 (10 - 2000 ng/ μ L)
Maximum Sample Concentration	2000 ng/ μ L
Concentration (CV) ^a	< 10%
Sizing Resolution ^b	\leq 10% difference in size
Sizing Precision RSD (CV)	< 2%
Separation Time per Sample	65 seconds
Sensitivity (LOD) ^c	0.2 ng/ μ L
Reagent Kit Primes	10
Chip Lifetime	HT: 400 samples 24: 400 samples
Samples per Calibrated Chip Prep	HT: up to 96 samples LT: up to 48 samples
HT Preps per Chip (96 samples)	4
LT Preps per Chip (48 samples)	8
Minimum Sample Volume	2 μ L

a. Typical results. Concentration CV may vary \pm 5% based on individual proteins.

b. Resolution is defined as the difference in migration times divided by the sum of the full width half max for two closely migrating peaks.

c. Based on internal standards.

Sample Conditions

Table 2. Sample Conditions

Buffers, Salts and Additives	Refer to “ Compatible Buffers, Salts and Additives ” on page 37 for compatibility with specific buffers, salts and additives. If your conditions are not listed, contact Revvity (see page 39) for more information on compatibility.
Particulates	Spin down sample plates prior to analysis. Filter all buffers with a 0.22 µm cellulose acetate filter.
Salt Concentration	Total salt concentration must not exceed 1M.

Storage Conditions

Chip Storage: Prior to use, store chips refrigerated at 2 - 8°C. After use, store chips at room temperature and use within 30 days.

Reagent Storage:

- Store the Protein Clear HR Dye solution (blue cap) and the ProteinEXact HR Ladder (yellow cap) at -20°C.
- Store all other reagents at 2 - 8°C.
- After preparation, store the Gel-Dye solution at 2 - 8°C.

CRITICAL:

The sample plate, chip, and all refrigerated reagents must equilibrate to room temperature (20 - 25°C) for at least 30 minutes before use. Protect the Protein Express Lower Marker from light when not in use.

Remove the Protein Clear HR Dye and the ProteinEXact HR Ladder from the padded shipping pack and allow to warm from -20°C to room temperature for 45 minutes. Protect the Protein Clear HR Dye from light.

To avoid multiple freeze/thaw cycles, aliquot 15 µL of the ProteinEXact HR Ladder into each of the provided empty ladder tubes. Store these aliquots, including the original tube, at -20°C.

Reagent Kit Contents

ProteinEXact HR Reagent Kit Contents P/N CLS150466

Table 3. Reagents

Reagent	Vial	Quantity
Protein Clear HR Dye solution	Blue 	1 vial, 0.13 mL
Protein Express Sample Buffer	White 	4 vials, 1.5 mL each
ProteinEXact Gel Matrix	Red 	3 vials, 1.7 mL each
ProteinEXact HR Ladder	Yellow 	2 vial, 0.08 mL each
Protein Express Lower Marker	Green 	2 vials, 0.5 mL each
Protein Express Wash Buffer	Purple 	5 vials, 1.8 mL each

Table 4. Consumable Items

Item	Supplier and Catalog Number	Quantity
Empty ladder tubes for aliquotes	Revvity, Cat. # CLS160196	6
Spin Filters	Costar [®] , Cat. # 8160	20
Detection Window Cleaning Cloth	VWR [™] , Cat. # 21912-046	1
Swab	ITW Texwipe [®] , Cat. # TX758B	3
Ladder tubes, 0.2 mL	(Not sold separately)	20
Buffer Tubes, 0.75 mL	(Not sold separately)	20

LabChips

Table 5. ProteinEXact HR LabChips

Item	Part Number
HT ProteinEXact HR LabChip (for use with GXII Touch HT)	CLS150337
24 ProteinEXact HR LabChip (for use with GXII Touch 24 or HT)	CLS150338

Safety and Usage

Safety Warnings and Precautions

CAUTION

We recommend that this product and components be handled only by those who have been trained in laboratory techniques and that it is used in accordance with the principles of good laboratory practice. As all chemicals should be considered potentially hazardous, it is advisable when handling chemical reagents to wear suitable protective clothing, such as laboratory overalls, safety glasses, and gloves. Care should be taken to avoid contact with skin or eyes. In case of contact with skin or eyes, wash immediately with water.

WARNING!



- *Dye solution contains DMSO. Avoid contact with skin and eyes.*
- *Dye solution contains SDS. Avoid inhalation and contact with skin and eyes.*
- *Wash Buffer and Sample Buffer contain LDS. Avoid inhalation and contact with skin and eyes.*
- *Gel Matrix contains Methyl Urea. Avoid contact with skin and eyes.*

Usage

The ProteinEXact HR Assay is for use with LabChip GXII Touch instruments. LabChip GXII Touch instruments are for research use only and not for use in diagnostic procedures.

Preparation Procedures

CRITICAL:

- *The sample plate, chip, and all refrigerated reagents must equilibrate to room temperature (20 - 25°C) for at least 30 minutes before use. Protect the Protein Express Lower Marker from light.*
- *Remove the ProteinEXact HR Ladder (yellow cap) from the padded shipping pack and allow to warm from -20°C to room temperature for 45 minutes. To avoid multiple freeze/thaw cycles, aliquot 15 µL of the ProteinEXact HR Ladder into each of the provided empty ladder tubes. Store these aliquots, including the original tube, at -20°C..*
- *Remove the Protein Clear HR Dye (blue cap) from the padded shipping pack and allow to warm from -20°C to room temperature for 45 minutes, protected from light.*
- *The assay requires exact and consistent adherence to the protocol as shown in this guide, or results may be compromised by increased variability.*
- *Fresh Milli-Q® water should be obtained the day of the assay.*
- *Adherence to the full vortex time is important for assay performance.*

Additional Items Required

- 0.6 mL centrifuge tubes and/or 96-well plates for denaturing protein samples.
- Means for heating samples to 70°C — 96-well PCR instrument or heating block.

Note: *Avoid using non-stick lab consumables. They may induce unexpected or erratic assay results caused by surface treatments leaching into dye or gel components.*

- 18 megohm, 0.22-µm filtered water (Milli-Q® or equivalent).
- 70% isopropanol solution in DI water.
- Reducing agents: BME (beta-mercaptoethanol), 1M DTT (dithiothreitol) or 100 mM TCEP.
- Non-reducing agents: 250 mM IAM (iodoacetamide)

Preparing the Gel-Dye Solution

Notes:

- *The dye is light sensitive. **Do not expose the Dye solution or Gel-Dye solution to light for any length of time.** Keep the prepared Gel-Dye solution in the dark.*
- *Store the Gel-Dye solution in the dark for 3 weeks at 2 - 8°C.*

For High Throughput Chip Preparation (up to 96 samples)

- 1 Vortex the thawed Protein Clear HR Dye solution at max speed for 20 seconds and quickly spin down before use.

Note: *Adherence to the full vortex time is important for assay performance.*

- 2 Using a reverse pipetting technique, transfer 520 μ L of ProteinEXact Gel Matrix (red cap ●) to the top “basket” of a provided spin filter.

Note: *Gel Matrix is extremely viscous. It is important to use a reverse pipetting technique, as described on [page 31](#), to accurately transfer the correct amount of gel to the spin filter. Incorrect ratios of gel to dye will cause inconsistent assay results.*

- 3 Add 20 μ L of Protein Clear HR Dye solution (blue cap ●) to the 520 μ L Gel Matrix in the spin filter. For best results, make fresh and use immediately.
- 4 After the Dye solution is added to the Gel Matrix, immediately cap and invert the spin filter 10 times to mix well and minimize dye concentrate interaction with filter material; then vortex tube, upside down, for 20 seconds until the gel and dye are well mixed.

Note: *Adherence to the full vortex time is important for assay performance.*

- 5 For Destain Solution, transfer 250 μ L of ProteinEXact Gel Matrix (red cap ●) to a second spin filter.
- 6 Spin the Gel-Dye solution and the Destain Solution at 9300 rcf for 8 minutes at RT. Ensure that the microcentrifuge is set to RT and the material has passed through the filter (spin longer if necessary), then discard the filter baskets and cap the tubes. Store in the dark until ready to use.

Note: *Do not exceed 9300 rcf when filtering Gel-Dye solution. Exceeding 9300 rcf will change the properties of the gel.*

For Low Throughput Chip Preparation (48 samples)

- 1 Vortex the thawed Dye solution at max speed for 20 seconds and quickly spin down before use.

Note: Adherence to the full vortex time is important for assay performance.

- 2 Using a reverse pipetting technique, transfer 280 μ L of ProteinEXact Gel Matrix (red cap ●) to the top “basket” of a provided spin filter.

Note: Gel Matrix is extremely viscous. It is important to use a reverse pipetting technique, as described on [page 31](#), to accurately transfer the correct amount of gel to the spin filter. Incorrect ratios of gel to dye will cause inconsistent assay results.

- 3 Add 10.7 μ L of Protein Clear HR Dye solution (blue cap ●) to the 280 μ L Gel Matrix in the spin filter. For best results, make fresh and use immediately.
- 4 After the Dye solution is added to the Gel Matrix, immediately cap and invert the spin filter 10 times to mix well and minimize dye concentrate interaction with filter material; then vortex tube, upside down, for 20 seconds until the gel and dye solution are well mixed.

Note: Adherence to the full vortex time is important for assay performance.

- 5 For Destain Solution, transfer 180 μ L of ProteinEXact Gel Matrix (red cap ●) to a second spin filter.
- 6 Spin the Gel-Dye solution and the Destain Solution at 9300 rcf for 8 minutes at room temperature. Ensure that the microcentrifuge is set to RT and the material has passed through the filter (spin longer if necessary), then discard the filter baskets and cap the tubes. Store in the dark until ready to use.

Note: Do not exceed 9300 rcf when filtering Gel-Dye solution. Exceeding 9300 rcf will change the properties of the gel.

Preparing the Ladder and Wash Buffer Tube

Note: The ProteinEXact HR Ladder should be kept frozen. It is recommended that you aliquot 15 μL of the ProteinEXact HR Ladder into each of the provided empty ladder tubes for individual use after thawing for the first time. Store the aliquots at $-20\text{ }^{\circ}\text{C}$.

- 1 Ensure the ProteinEXact HR Ladder (yellow cap ●) has been warmed to room temperature, then vortex gently for 10 seconds. Briefly spin the ladder vial. Ensure no precipitate is visible in the solution. If precipitate is present, let the vial sit at room temperature for a little longer then repeat the gentle vortex and spin.

Note: Adherence to the full vortex time is important for assay performance.

- 2 Pipette 15 μL of ProteinEXact HR Ladder into a supplied Ladder Tube.
- 3 Add 150 μL of water (Milli-Q[®] or equivalent) to the Ladder Tube and mix thoroughly by pipetting up and down. Ensure that no bubbles are in the ladder tube.
- 4 Insert the Ladder Tube into the ladder slot on the LabChip GXII Touch instrument.
- 5 Transfer 750 μL of Protein Express Wash Buffer (purple cap ●) to the 0.75 mL Buffer Tube provided with the reagent kit. Ensure there are no air bubbles in the Buffer Tube.
- 6 Insert the Buffer Tube into the buffer slot on the LabChip GXII Touch instrument.

Note: Replace the Buffer Tube with a freshly prepared tube every 8 hours when the chip and instrument are in use.

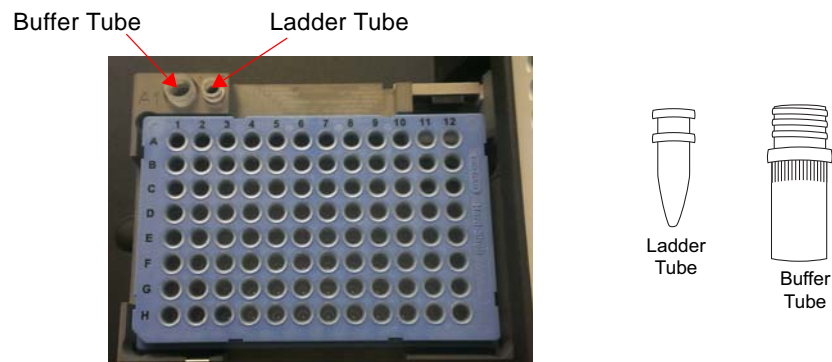


Figure 1. Buffer Tube and Ladder Tube in the GXII Touch

Preparing the Chip

- 1 Allow the chip to equilibrate to room temperature (20 - 25°C) for at least 30 minutes before use.
- 2 Use a pipette tip attached to a vacuum line to thoroughly aspirate all fluid from the chip wells (see [Figure 2](#)). For details on how to set up a vacuum line, see [page 38](#).
- 3 Rinse and completely aspirate each active chip well (1, 2, 3, 4, 7, 8, 9, and 10) two times with water (Milli-Q[®] or equivalent). Do not allow active wells to remain dry.
- 4 If any water spills onto the top or bottom of the chip surfaces during rinsing, aspirate using the vacuum line. DO NOT run the tip over the central region of the detection window. Use the provided Detection Window Cleaning Cloth to clean the chip detection window.

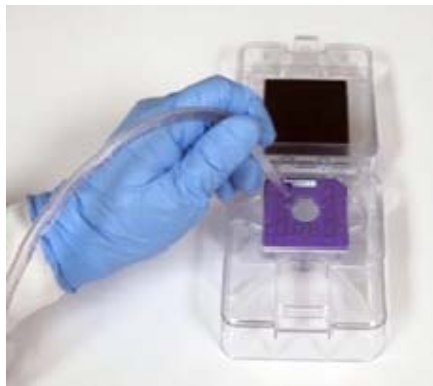


Figure 2. Using a vacuum to aspirate the chip wells is more effective than using a pipette. See [page 38](#) for details.

- 5 Using a reverse pipetting technique, add Gel-Dye solution from the spin filter tube to chip wells 3, 7, 8, and 10 with volumes shown in [Figure 3](#) (High-throughput) or [Figure 4](#) (Low-throughput).

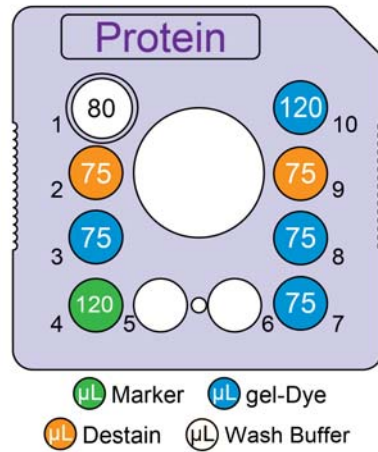


Figure 3. Reagent placement for High-throughput (up to 96 samples)

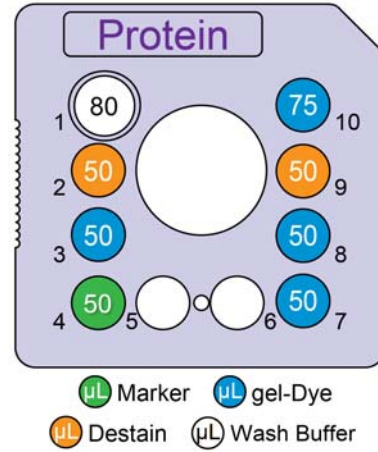


Figure 4. Reagent placement for Low-throughput (up to 48 samples)

- 6 Using a reverse pipetting technique, add Destain Solution from the spin filter tube to chip wells 2 and 9 with volumes shown in [Figure 3](#) (High-throughput) or [Figure 4](#) (Low-throughput).
- 7 Using a reverse pipetting technique, add Protein Express Lower Marker (green cap ●) to chip well 4 with volumes shown in [Figure 3](#) (High-throughput) or [Figure 4](#) (Low-throughput). Make sure the marker volume is pipetted accurately. If there is not enough marker in chip well 4, the marker will deplete and will not be added to subsequent samples on-chip. Data collected without marker peaks cannot be analyzed by the software.
- 8 Using a reverse pipetting technique, add 80 µL Wash Buffer (purple cap ●) to chip well 1 (waste well) for both High-throughput and Low-throughput as shown in [Figure 3](#) and [Figure 4](#).
- 9 Remove any liquid from the chip surface and the rims of the wells using the vacuum line.

Inserting a Chip into the LabChip GXII Touch

- 1 Check that the sample plate, Buffer Tube, and Ladder Tube are placed on the instrument properly.
- 2 Touch the *Load Plate* button on the *Home* screen (Figure 5) to retract the sample plate and move the sipper to the Buffer Tube.
- 3 Remove the chip from the storage container and inspect the detection window. Clean BOTH sides of the detection window with the Revvity-supplied Detection Window Cleaning Cloth dampened with a 70% isopropanol solution in DI water.
- 4 Touch the *Unload Chip* button on the *Home* screen.

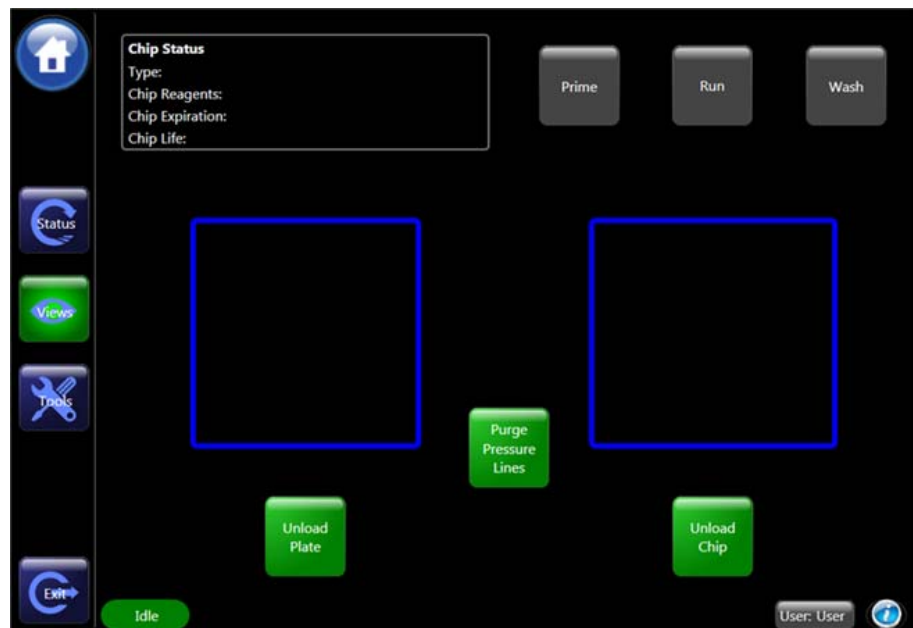


Figure 5. Home screen

- 5 Insert the chip into the LabChip GXII Touch instrument (Figure 6) and close the chip door securely.

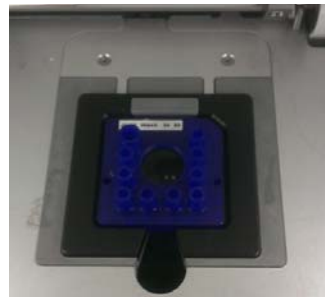


Figure 6. Chip in the LabChip GXII Touch instrument

Priming and Warming

Chips can be primed independently from running assays on the LabChip GXII Touch instrument. As a result, samples can be prepared during the priming process.

To prime the chip while preparing samples:

- 1 Place buffer tube, ladder tube, and chip in the instrument.
- 2 Touch the *Prime* button on the *Home* screen (Figure 7).

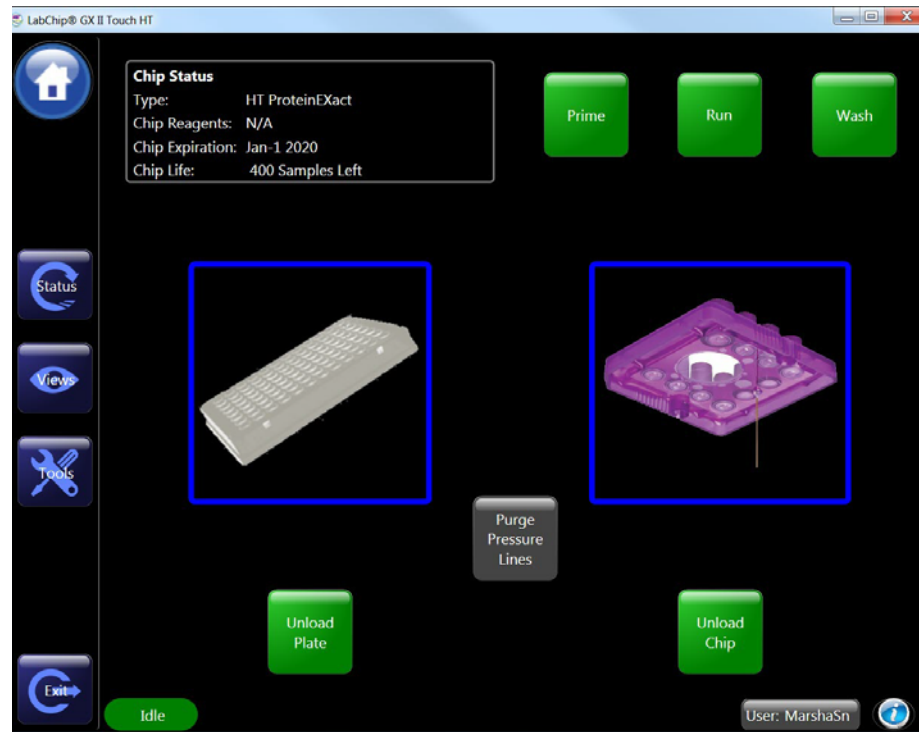


Figure 7. *Prime* button on *Home* screen

The *Prime* screen opens (Figure 8).

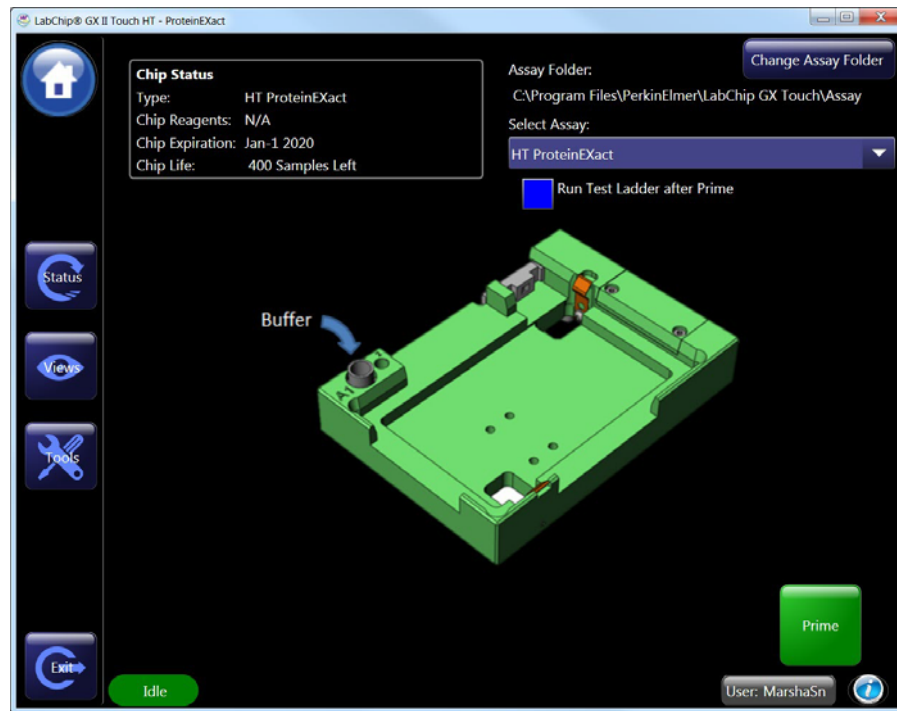


Figure 8. Prime button on the Prime screen

- 3 If desired, select the *Run Test Ladder after Prime* check box to run one ladder after the prime is complete.
- 4 Touch the *Prime* button to begin the prime and warming process.

When starting the run, the *Setup Run* tab will display a *Skip Warm* check box and a *Skip Prime* check box during the run (see [Figure 11 on page 18](#)) so the run can be started without repeating the priming and warming steps.

Note:

- *Do not keep the chip door open for any length of time after the chip is installed.*

Preparing the Protein Samples

Note: The following is a general protocol for antibody sample preparation. Optimization of the type or concentration of reducing agent and stabilizing agent and/or optimization of denaturing conditions may be necessary depending on the specific molecules to be analyzed. Use of a hardshell 96-well PCR plate and a thermal cycler is recommended for efficient sample preparation.

- 1 Prepare Reducing and/or Non-Reducing Sample buffer.
- 2 The sample buffer calculation is described as follows:
 - Transfer 700 μL of Protein Express Sample Buffer (white cap \bigcirc) into a microfuge tube.
 - For Reducing Sample Buffer, add 24.5 μL of BME or 1M DTT.
 - For Non-Reducing Sample Buffer, add 24.5 μL of 250 mM IAM.
- 3 For each sample to be analyzed, pipette 18 μL of Reducing or Non-Reducing Sample Buffer into a well in a 96-well PCR plate.
- 4 Add 2.5 μL of sample to each prepared well.
- 5 When finished, seal the plate and denature samples at 70°C for 10 min, then cool to room temperature. Optimum denaturing conditions may vary by sample type.
- 6 Add 35 μL of water (Milli-Q[®] or equivalent) to each sample and mix thoroughly by pipetting up and down.
- 7 Spin the sample plate at 3000 rpm for 2-3 minutes to eliminate bubbles and move the fluid to the bottom of the wells.
- 8 Place the sample plate onto the instrument's plate holder.

Running the Assay

- 1 Insert the sample plate, buffer, and ladder tubes.
- 2 Touch the *Run* button on the *Home* screen.
- 3 On the *Select Wells* tab (Figure 9), select the appropriate plate type, well pattern, and whether to read wells in columns or rows. Select number of times each well is sampled under *Adv. Settings*. Touch the green arrow.

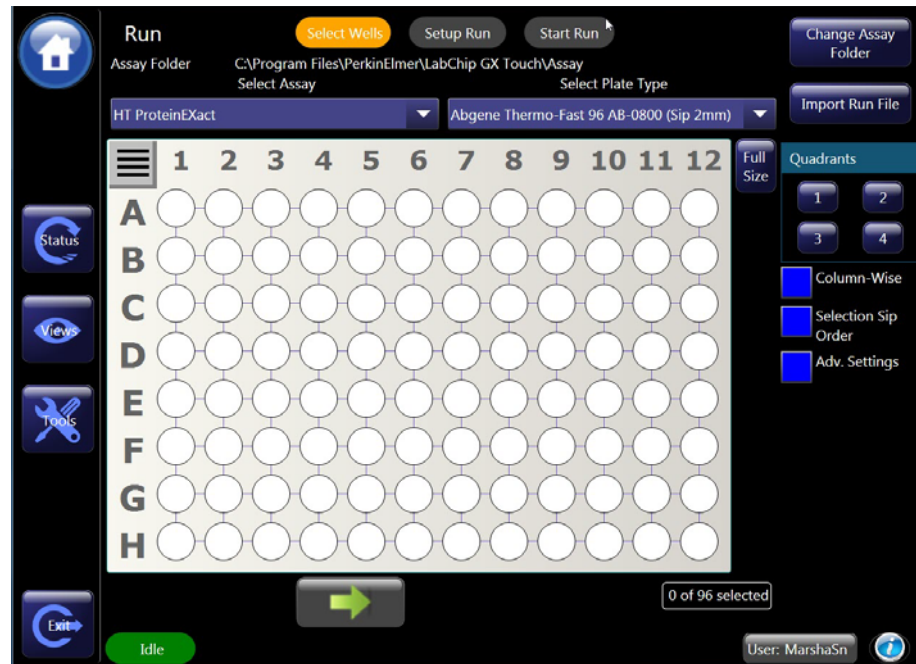


Figure 9. *Select Wells* Tab

- 4 If the chip has *not* yet been primed, as described in “[Priming and Warming](#)” on page 14, the *Setup Run* tab in [Figure 10](#) displays.



Figure 10. *Setup Run* tab when priming and warming takes place during run

If the chip was primed and warmed before the run, the *Skip Warm* and *Skip Prime* check boxes display as shown in [Figure 11](#).



Figure 11. *Setup Run* tab when priming and warming takes place before the run

- 5 Select the operator name, the option to read barcode, the destination of the file, the inclusion of sample names, expected peaks, excluded peaks, and the desired filename convention.
- 6 If the chip was primed and warmed before the run, select the *Skip Prime* and *Skip Warm* check boxes to skip priming at the beginning of the run.

- 7 If desired, select *Auto Export* to export results tables automatically, *Auto Print* to print the results table to a PDF, and *Defer Export to Plate Completion* to export the results at the end of the run instead of as each well is completed.
- 8 Touch the green arrow. The *Start Run* tab displays (Figure 12).



Figure 12. *Start Run* Tab

- 9 Type the total ladder concentration (in mg/mL) in the *Total Ladder Concentration* text box. The Ladder Concentration is specified on the ProteinEXact HR Ladder vial label as shown in Figure 13. The ladder concentration should be within the range of 0.499 mg/mL +/- 15%.

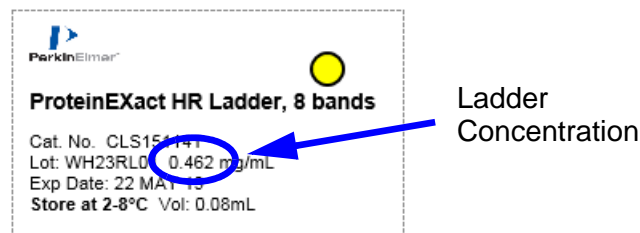


Figure 13. Ladder Concentration on ProteinEXact HR Ladder Vial Label

- 10 Touch the *Start* button to begin the run.

Cleaning and Storing the Chip

After use, the chip must be cleaned and stored in the chip container.

- 1 Place the chip into the chip storage container. The sipper must be submerged in the fluid reservoir.
- 2 Remove the reagents from each well of the chip using vacuum.
- 3 Rinse and completely aspirate each active well (1, 2, 3, 4, 7, 8, 9, and 10) two times with water (Milli-Q[®] or equivalent).
- 4 Add 120 μ L water (Milli-Q[®] or equivalent) to the active wells.
- 5 Cover the wells with Parafilm[®] to prevent evaporation and store the chip at room temperature. Allowing chip wells to dry may lead to changes in chip performance. The chip must be used to its lifetime (to the total number of 400 samples) within 30 days of analyzing the first plate of samples.

Chip Cartridge Cleaning

1 Daily Cleaning

- a Inspect the inside of the chip cartridge and O-rings for debris.
- b Touch the *Purge Pressure Lines* button on the Home screen (see [Figure 5 on page 13](#)).
- c Use the provided lint-free swab dampened with water (Milli-Q[®] or equivalent) to clean the electrodes and the O-rings using a circular motion. If the O-rings stick to the chip or if a pressure leak is detected, perform the more extensive Monthly cleaning procedure.

2 Monthly Cleaning

- a To reduce pressure leaks at the chip interface, clean the O-rings frequently. Remove the O-rings from the top plate of the chip interface on the LabChip GXII Touch instrument. Soak O-rings in water (Milli-Q[®] or equivalent) for several minutes. Clean the O-ring faces by rubbing between two fingers. Wear gloves.
- b To reduce the occurrence of current leaks, clean the chip interface frequently. Clean the top plate of the chip interface using the provided lint free swab dampened with water (Milli-Q[®] or equivalent).
- c Allow the O-rings and chip interface to air dry. Reinsert the O-rings into the chip cartridge.

Results

ProteinEXact HR Ladder Result

The electropherogram of a typical ProteinEXact HR ladder is shown in [Figure 14](#). Peaks to the right of the lower marker and system peaks in order of increasing migration time correspond to proteins of increasing size i.e., 6.5 kDa, 15.9 kDa, 20.4 kDa, 28.9 kDa, 48.4 kDa, 68.4 kDa, 119.2 kDa, and 250 kDa, respectively.

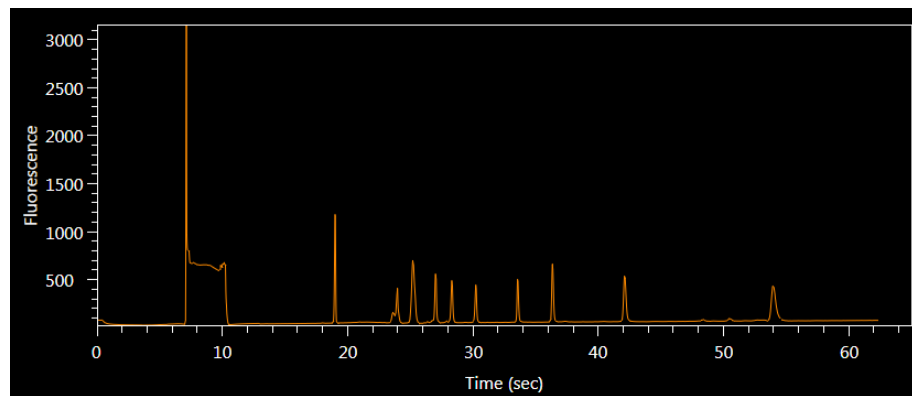


Figure 14. ProteinEXact HR ladder electropherogram

Troubleshooting

Note: Some of the data examples in this section were generated with assays other than the assay described in this user guide.

Symptom: Failed calibration.

Possible causes:

- 1 Standard area measurement is outside assay validation range. The instrument gain factor has shifted.
- 2 Expected standard arrival time is outside assay validation range. Gel-dye mixing ratios are different from nominal or a poor chip prime.

What to do:

- 1 Calibrate the optics as described in the *LabChip GX Touch/LabChip GXII Touch User Manual*.

Note: Software versions 1.9 and later select the minimum or maximum laser power if standard area measurement is outside of assay validation range and will not display as a calibration failure.

- 2 Re-prepare the chip, making sure the reagents are fresh, and retry calibration by performing another run.

Symptom: No ladder or sample peaks but marker peaks detected.

Note: The lower marker peak height will most likely be greater than normal height.

Possible causes:

- 1 Air bubble or clog in sipper introduced during chip priming.

What to do:

- 1 Reprime the chip. See [“Repriming Chips” on page 33](#) for instructions on how to reprime the chip.
- 2 Perform a sipper unclogging procedure. See [“Removing Sipper Clogs” on page 36](#) for instructions.

Symptom: Missing sample, ladder *and* marker peaks.**Possible causes:**

- 1 Clog in sipper or marker channel of chip.

What to do:

- 1 Reprime the chip. See [“Repriming Chips” on page 33](#) for instructions on how to reprime the chip.

Symptom: Ladder detected but no sample peaks.**Possible causes:**

- 1 The sipper is not reaching the sample due to low sample volume in the well of the plate.
- 2 If the missing sample peaks occurred only in a few wells of the plate, check those wells for air bubbles.
- 3 The sipper is not reaching the sample due to an incorrect capillary height setting or incorrect plate definition.
- 4 If the plate has been uncovered for some time, sample evaporation might have occurred.
- 5 Debris from the sample or sample prep is clogging the sipper.

What to do:

- 1 Add more sample to the well.
- 2 Manually insert a larger volume pipette tip (~100 μ L) into the sample well and dislodge the bubble. Rerun these sample wells.
- 3 Check the plate definitions.
- 4 Check the sample wells, especially around the edge of the plate where evaporation is fastest, and make a fresh plate if volumes are low.
- 5 If you suspect there may be debris in your samples, spin the sample plate down in a centrifuge (e.g., 3000 rpm for 5 minutes). Unclog the sipper by repriming the chip. See [“Repriming Chips” on page 33](#) for instructions on how to reprime the chip.

Symptom: No ladder peaks but sample peaks and marker peaks are present.**Possible causes:**

- 1 Low or no ladder volume in the Ladder Tube.

What to do:

- 1 Add more ladder to the Ladder Tube and restart the run. Recommended standard ladder volume is 120 μ L (minimum volume is 100 μ L).

Symptom: No marker peaks but sample peaks are present.**Possible causes:**

- 1 No marker added to chip well 4.
- 2 If there is marker solution in chip well 4, the problem may be due to a marker channel clog.

What to do:

- 1 This may be due to not filling the marker well or the chip remaining idle on the instrument for an extended period of time. Add or replenish the Lower Marker in the chip using the following procedure:
 - Touch the *Unload Chip* button on the Home screen to open the chip door.
 - Return the chip to the chip container ensuring the sipper is immersed in fluid.
 - Thoroughly aspirate all fluid from chip well 4 using a vacuum line.
 - Rinse and completely aspirate chip well 4 twice with water (Milli-Q[®] or equivalent).
 - Add Protein Express Lower Marker (green cap ●) to chip well 4.
 - Insert the chip back into the instrument.
 - Restart the run.
- 2 Perform a marker channel unclogging procedure by repriming the chip. See [“Repriming Chips” on page 33](#) for instructions on how to reprime the chip.

Symptom: Ladder traces show up in the lanes following the ladders (delayed sip).

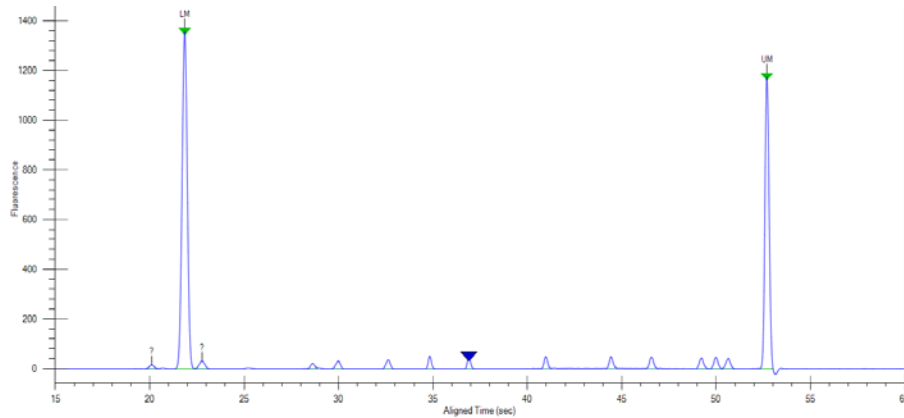


Figure 15. Small ladder peaks in sample well caused by delayed sip

Possible causes:

- 1 Separation channel overloaded with sample.
- 2 Partial clog in the separation channel.

What to do:

- 1 Lower the starting sample concentration.
- 2 Reprime the chip. See [“Repriming Chips” on page 33](#) for instructions on how to reprime the chip.

Symptom: Unexpected sharp peaks.

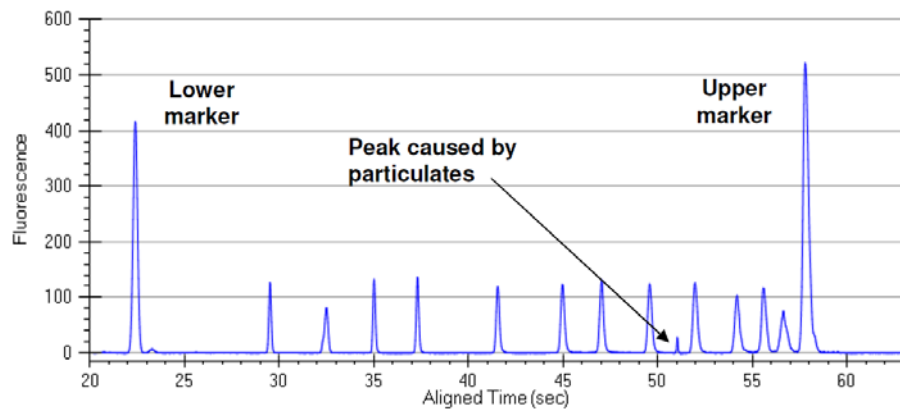


Figure 16. Unexpected sharp peak

Possible causes:

- Dust or other particulates introduced through sample or reagents.

What to do:

- 1 Do one or all of the following:
 - Replace the 18 megohm, 0.22- μm filtered water (Milli-Q[®] or equivalent) water used for chip preparation.
 - Replace the buffer used for sample and reagent preparation.
 - Use a 0.22-micron filter for all water and buffers used for chip, sample, and reagent preparation.
 - Spin down sample plate to pellet any particulates.

Symptom: Humps in several electropherograms which do not correspond to sample data.

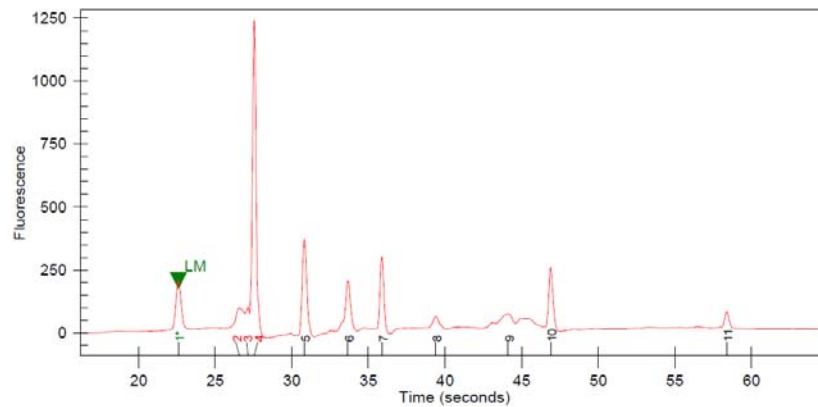


Figure 17. Humps in several electropherograms

Possible causes:

- 1 Electrode 7 is dirty and has contaminated the Gel-Dye solution in well 7.
- 2 High concentrations of detergent in the sample buffer can sometimes cause humps in the electropherogram.

What to do:

- 1 Before restarting the run, clean electrode 7. Remove the chip and follow the electrode cleaning procedure. We recommend using the provided swab and isopropanol to manually clean electrode 7.
- 2 Lower the detergent concentration in the sample (see [“Compatible Buffers, Salts and Additives”](#) on page 37).

Symptom: Peaks migrating much faster or slower than expected.

Note: Some migration time variance between chips or within a plate is considered normal chip performance. All chips are QC tested at Revvity prior to shipment.

Normal migration time windows for the markers are:

- ProteinEXact HR Assay Lower Marker: 18-19 seconds
- 120 kD Ladder Protein on the first plate: 41-45 seconds

Possible causes:

- 1 Incorrect Gel-Dye ratio. Migration time is sensitive to dye concentration and peaks will migrate too fast or too slow if the dye concentration in the gel is too low or too high, respectively.

Note: Excess dye in the separation channel will slow down migration, and less dye in the separation channel will make peaks migrate faster.

- 2 Particulates from the samples may be clogging the separation channel (this will slow down migration).
- 3 Gel-Dye solution was not primed properly into the chip.

What to do:

- 1 Prepare fresh Gel-Dye solution. Wash and reprime the chip with the new Gel-Dye solution. See [“Washing Chips” on page 34](#) for instructions on how to wash and reprime the chip.
- 2 If fast or slow migration is observed repeatedly on a new chip, contact technical support to arrange return of the chip to Revvity. Please send a data file showing the failure along with the return request.
- 3 Minimize the loading of particulates in the sample by performing a centrifuge spin of the sample plate (e.g. 3000 rpm for 5 minutes) and/or ensuring the Sip 4 mm plate type is selected in the Select Wells screen before starting a new run. The debris may be flushed out of the chip by washing and re-priming the chip. See [“Washing Chips” on page 34](#) for instructions on how to wash and reprime the chip.
- 4 Check the O-rings on the top surface of the chip interface and clean if necessary.

Symptom: High baseline fluorescence (e.g., greater than 1000 counts).**Possible causes:**

- 1 The destain wells (2 and 9) do not contain Destain solution (Gel Matrix with no dye).
- 2 The destain wells (2 and 9) may have been contaminated with dye either because the well was improperly flushed after priming or because dye solution was mistakenly pipetted into the well.

What to do:

- 1 Prepare a fresh Destain solution. Wash and reprime the chip with the new Destain solution. See [“Washing Chips” on page 34](#) for instructions on how to wash and reprime the chip.

Symptom: Lower than expected signal for ladders and samples.**Possible causes:**

- 1 Improper SDS concentration in Gel-Dye solution.

What to do:

- 1 Ensure that Dye solution is completely thawed and mixed. Prepare fresh Gel-Dye solution. Wash and reprime the chip with the new Gel-Dye solution. See [“Washing Chips” on page 34](#) for instructions on how to wash and reprime the chip.

LabChip Kit Essential Practices

To ensure proper assay performance, please follow the important handling practices described below. Failure to observe these guidelines may void the LabChip Kit product warranty.¹

Note: *It is important to keep particulates out of the chip wells, channels, and capillary. Many of the following guidelines are designed to keep the chips particulate-free.*

For assay and instrument troubleshooting, refer to the Software Help file or call Revvity Technical Support at 1-800-762-4000.

General

- The assay requires exact and consistent adherence to the protocol as shown in this guide, or results may be compromised by increased variability.
- Warm the sample plate, chip, and reagents to room temperature (20 - 25°C) as described in [“Preparation Procedures” on page 7](#)
- Clean the chip interface O-rings weekly and the electrodes daily. Refer to the Instrument Users Guide Maintenance and Service section for procedures.
- Avoid use of powdered gloves. Use only non-powdered gloves when handling chips, reagents, sample plates, and when cleaning the instrument electrodes and electrode block.
- Calibrate laboratory pipettes regularly to ensure proper reagent dispensing.
- Only the Revvity-supplied Detection Window Cleaning cloth can be used on the chip to clean the detection window.
- Water used for chip preparation procedures must be 18 megohm, 0.22-µm filtered water (Milli-Q® or equivalent).
- Using the “Reverse Pipetting Technique” (see [page 31](#)) will help avoid introducing bubbles into the chip when pipetting the gel.

1. Revvity, Inc. warrants that the LabChip Kit meets specification at the time of shipment, and is free from defects in material and workmanship. LabChip Kits are warranted for 60 days from the date of shipment. All claims under this warranty must be made within thirty days of the discovery of the defect.

Reverse Pipetting Technique

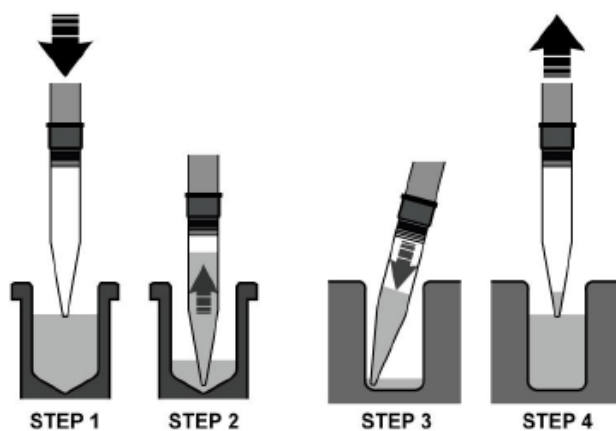


Figure 18. Reverse pipetting

- 1** Depress the pipette plunger to the first stop, then continue to depress the plunger past the first stop, but not all the way to the second stop.
- 2** Aspirate the selected volume plus an excess amount from the tube. Leave the pipette tip in the solution for 5-10 seconds to ensure that the aspiration is complete.
- 3** Dispense the selected volume into the corner of the well by depressing plunger to the first stop. Hold the pipette for 10 seconds with the plunger at the first stop to ensure that the dispense is complete.
- 4** Withdraw the pipette from the well.
- 5** Retain the excess volume in the pipette tip if performing an additional Reverse Pipetting transfer, or return the excess volume to the source container for later use.

Reagents

- Store reagents as specified in [“Storage Conditions” on page 4](#).
- Warm the sample plate, chip, and reagents to room temperature (20 - 25°C) as described in [“Preparation Procedures” on page 7](#)
- Protect the Dye, Gel-Dye solution, and Protein Express Lower Marker from light. Store in dark when not in use.
- The Gel-Dye solution expires 3 weeks after preparation.
- For optimal performance, use one reagent kit per chip. The ProteinEXact HR Reagent Kit contains the reagents to run 400 samples, equivalent to four chip preparations in the high-throughput mode or eight chip preparations in low-throughput mode, whichever comes first.

Chips

General Guidelines

- Store chips as specified in [“Storage Conditions” on page 4](#).
- Warm the chip to room temperature (20 - 25°C) as described in [“Preparation Procedures” on page 7](#) before use.
- After use, store chips at room temperature and use within 30 days
- Do not allow the liquid in the chip container to freeze, as this may lead to poor chip performance. Do not submerge the chip in any solution.
- The entire chip surface must be thoroughly dry before use.
- The sipper must be kept immersed in fluid at all times and should not be exposed to an open environment for long periods of time.
- Use care in chip handling to prevent sipper damage. Damage to the sipper can result in inconsistent sampling.
- Avoid exposing the chip to dust by keeping the chip in a closed environment such as in the chip container or in the instrument before and after chip preparation.
- Chips can be prepared and left in the instrument for extended periods of time so that samples can be run as needed throughout the day. Revvity recommends the chip be re-prepared after it has been idle for 8 hours, but the chip can be used continually over an 8-hour work day as long as the maximum recommended idle time of 8 hours and total chip lifetime of 400 samples are not exceeded.

Repriming Chips

- 1 Touch the *Unload Chip* button on the *Home* screen to open the instrument door.
- 2 Place the chip into the instrument.
- 3 Close the chip door securely and choose the corresponding assay.
- 4 Touch the *Prime* button on the *Home* screen and then touch the *Prime* button on the *Prime* screen.

OR

Click the *Start* button on the *Home* screen and then clear the *Skip Prime* and *Skip Warm* check boxes on the *Setup Run* tab.

Washing Chips

Important Note: Wash chips only with water (Milli-Q® or equivalent). Use of any other reagents (including Wash Buffer) is likely to cause even more artifacts in subsequent data.

Notes: Some protein samples may have components which produce data with extra peaks, spikes or other artifacts. When these artifacts are present, washing chips on the LabChip GXII Touch immediately before the next use can often restore data quality.

Chips should only be washed on the LabChip GXII Touch immediately before they are prepared with fresh reagents and primed on the instrument. Chips should not be washed and left with water in the chip channels for any extended period of time.

For most protein samples, the only chip cleaning protocol that is required is to rinse and aspirate the active wells twice with water (Milli-Q® or equivalent), and store the chip with 120 µL of water in each active well.

To wash a chip immediately before running an assay:

- 1 Thoroughly aspirate all fluid from the chip wells using a vacuum line.
- 2 Rinse and completely aspirate each active well (1, 2, 3, 4, 7, 8, 9, and 10) two times with water (Milli-Q® or equivalent). Do not allow active wells to remain dry.
- 3 Add 120 µL of water (Milli-Q® or equivalent) to each active well (1, 2, 3, 4, 7, 8, 9, and 10).
- 4 Touch the *Unload Chip* button on the *Home* screen and place the chip into the instrument.
- 5 Close the chip door securely.
- 6 Add 750 µL of water (Milli-Q® or equivalent) into the Buffer Tube. Install into the instrument.
- 7 Touch the *Wash* button on the *Home* screen (Figure 5 on page 13). The wash cycle begins (Figure 19).

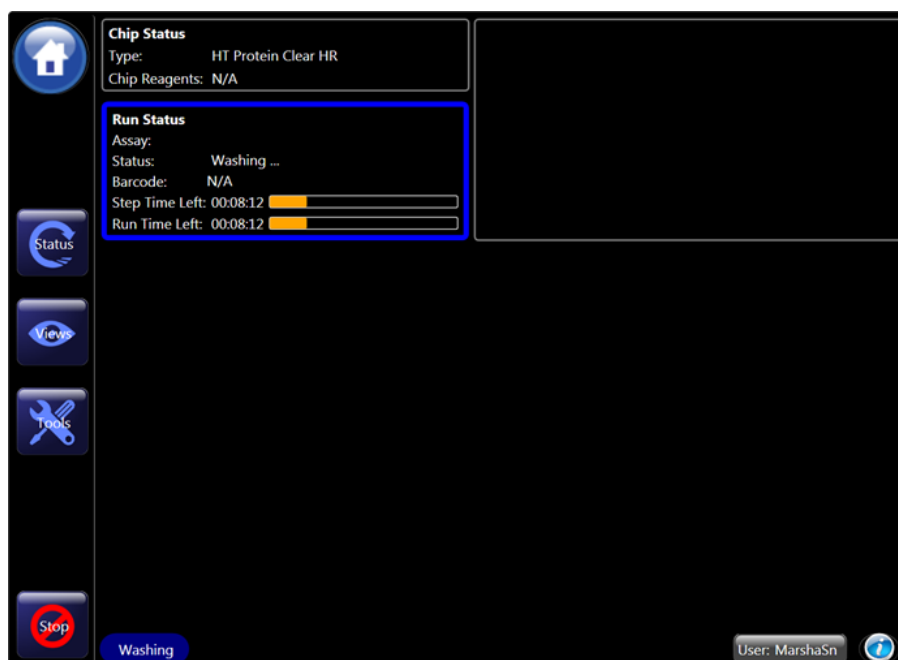


Figure 19. Wash screen

- 8 After completion of the wash cycle, open the chip cartridge and return the chip to the chip container, ensuring the sipper is immersed in liquid.
- 9 Thoroughly aspirate all liquid from the chip wells using a vacuum line.
- 10 Replace liquid in the wells with freshly made reagents as described in [“Preparing the Chip” on page 11](#). Do not let wells remain dry.
- 11 Transfer 750 μL of Wash Buffer (purple cap ●) into the supplied 0.75 mL Buffer Tube. Install into the instrument.
- 12 Install the Ladder Tube and sample plate in the instrument.
- 13 Run the assay.

Removing Sipper Clogs

If air bubbles are not dislodged after a reprime:

- Remove the chip from the instrument
- Rinse and completely aspirate well 1 (the waste well) twice with water (Milli-Q[®] or equivalent).
- Use reverse pipetting (described on [page 31](#)) to add 100 μL water (Milli-Q[®] or equivalent) to well 1.
- Suction the sipper with a vacuum line, as shown in [Figure 20](#), until droplets of fluid flow out from the sipper. When suctioning the sipper, be careful not to bend or break the sipper. To facilitate this, cut the end of the pipette tip attached to the vacuum line to widen the mouth.
- Aspirate the water from well 1, reinstall the chip into the instrument, and restart the run.

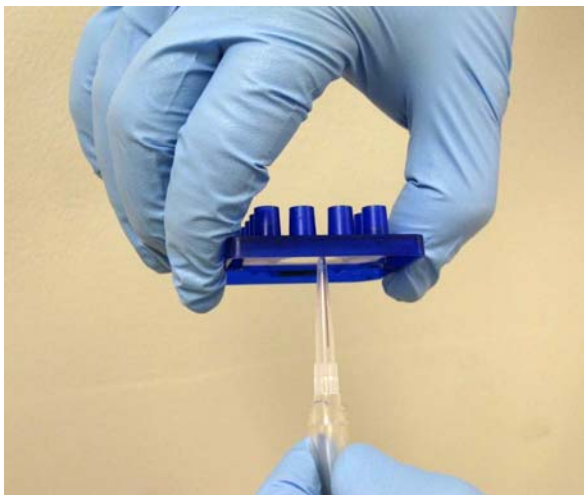


Figure 20. Removing an air bubble or clog by suctioning the sipper with a vacuum line

Samples

- Prepared sample plates should be free of gas bubbles and particulate debris, both of which may inhibit sipper flow.
- Sample plates containing gas bubbles and/or particulate debris should be spun down at 3000 rpm (1250 rcf) prior to analysis.
- Up to one 96-well plate (96 samples) can be run with a single HT chip preparation when running the ProteinEXact assay on the GXII Touch HT instrument.
- Up to 48 samples can be run with a single LT chip prep.

Compatible Buffers, Salts and Additives

Table 6. Compatible Buffers, Salts and Additives

Buffer & Salts	Concentration Limit	Additives	Concentration Limit
Tris Chloride	250 mM	Octyl Glucoside	2.5%
Tris Glycine	250 mM	Pluronic F68	0.1%
HEPES	500 mM	Sarcosyl	10%
PBS	8 X	CHAPS	0.5%
Sodium Citrate	150 mM	Tween 20	0.8%
Sodium Phosphate	250 mM	Triton X-100	0.6%
Sodium Acetate	600 mM	SDS	2%
Sodium Chloride	1000 mM	Zwittergent 3-14	0.4%
Sodium Azide	6%	PEG 3350	1%
Sodium Hydroxide	500 mM	Glycerol	30%
Potassium Chloride	900 mM	Urea	8 M
Ammonium Bicarbonate	1000 mM	Sucrose	1 M
Magnesium Chloride	300 mM	DMSO	25%
Imidazole	900 mM	EDTA	100 mM
PhosphoSafe		Ethanol	50%
BugBuster	2.5 X		
BPER			
POP Culture			
Insect POP Culture			

Table 7. Incompatible Buffers, Salts and Additives

Buffer & Salts	Concentration Limit	Additives	Concentration Limit
RIPA	All	None	N/A

Chip Well Aspiration Using a Vacuum

Aspirating with a pipette can leave used reagents in the chip wells. For this reason, Revvity recommends vacuuming the wells instead. This can be accomplished by attaching a permanent pipette tip to a house vacuum line with trap (Figure 21). To avoid contamination, use a new disposable pipette tip over the permanent tip for each chip aspirated (Figure 22).

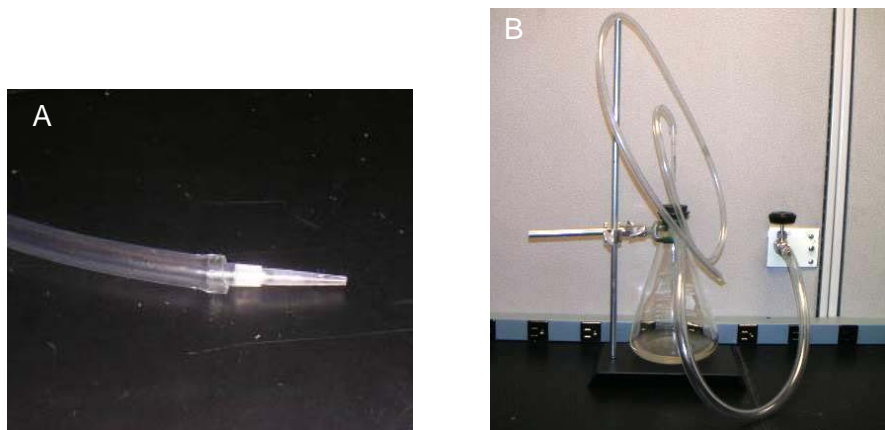


Figure 21. A: Permanent pipette tip attached to a house vacuum line; B: vacuum line with trap



Figure 22. Replacing the disposable pipette tip

Customer Technical Support

Revvity, Inc.
68 Elm Street
Hopkinton, MA 01748-1668 USA

Revvity Technical Support
Phone (USA Toll Free): 800-762-4000
Phone (Worldwide): +1 203-925-4602
Fax: +1 203-925-4602
Email: L3LabChip@Revvity.com
Internet: www.Revvity.com

For additional assay and instrument troubleshooting, refer to the Software Help file.

Licenses and Rights of Use

Label Licenses

This product is provided under an intellectual property license from Life Technologies Corporation. The purchase of this product conveys to the buyer the non-transferable right to use the data for internal research and training purposes solely in combination with Revvity, Inc. instrumentation, and not to generate revenue, which may include, but is not limited to use of the product: (1) in manufacturing or in quality assurance or quality control of a commercial product; (2) to provide a service, information, or data for a fee or other consideration; (3) for therapeutic or prophylactic purposes; (4) for diagnostic use; and (5) for resale, whether or not such items are resold for use in research. For information on purchasing a license to this product for any other purposes, contact Life Technologies Corporation, 5791 Van Allen Way, Carlsbad, CA 92008 USA or outlicensing@lifetech.com.

Rights of Use

The chip and reagents supplied with this kit are sold with limited rights of use. The chip may only be used with the specific quantity of reagents supplied with this kit. The purchaser has no right or license to refurbish, reuse, remanufacture, or otherwise use the chip with any other reagents than those specifically supplied in this kit. For more information on the terms and conditions of use of these chips and reagents, please read your LabChip GX Touch User Guide and refer to the applicable label license. The reagent kit contains materials that are provided under a license from Life Technologies Corporation, for research use only.

Trademarks

Revvity is a registered trademark of Revvity, Inc. All other trademarks are the property of their respective owners.

Copyright

This manual is published by Revvity, Inc., 68 Elm Street, Hopkinton, MA 01748 USA. Copyright 2018-2023 Revvity, Inc. All rights reserved.



Revvity, Inc.
68 Elm Street
Hopkinton, Massachusetts 01748 U.S.A.
TEL 508-435-9500
FAX 508-435-3439
<http://www.Revvity.com>

