



HTRF Insulin Ultra-Sensitive Detection Kit

Part # 62IN2PEG & 62IN2PEH

Test size#: 500 tests (62IN2PEG) and 10,000 tests (62IN2PEH) - assay volume: 20 μ L

Revision: #06 of October 2025

Store at: 2-8°C (62IN2PEG); 2-8°C (62IN2PEH)

For research use only. Not for use in diagnostic procedures.

ASSAY PRINCIPLE

This kit is intended for the simple and rapid quantification of Insulin in slightly concentrated pancreatic beta cell supernatants/lysates, islet supernatants/lysates and pancreas perfusates and offers a fast alternative to ELISA. The assay is compatible with human, mouse, rat, porcine and bovine species, but is not designed for use with plasma or serum samples.

The detection principle of this kit is based on HTRF® technology (Homogeneous Time-Resolved Fluorescence). As shown in Figure 1, Insulin is detected in a sandwich assay by using anti Insulin antibody labeled with Europium cryptate (donor), and anti Insulin antibody labeled with XL665 (acceptor).

When the dyes are in close proximity, the excitation of the donor with a light source (laser or flash lamp) triggers a Fluorescence Resonance Energy Transfer (FRET) towards the acceptor, which in turn fluoresces at a specific wavelength (665 nm). Signal intensity is proportional to the number of antigen-antibody complexes formed and therefore to the Insulin concentration.

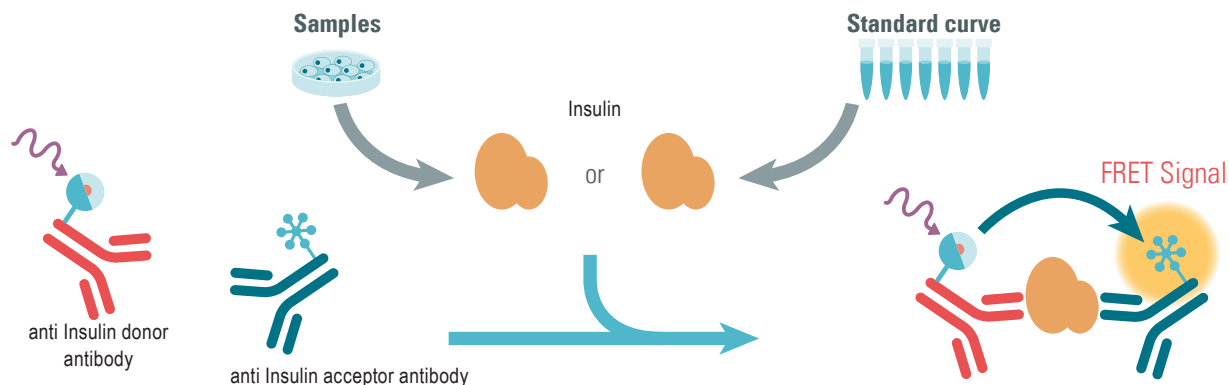
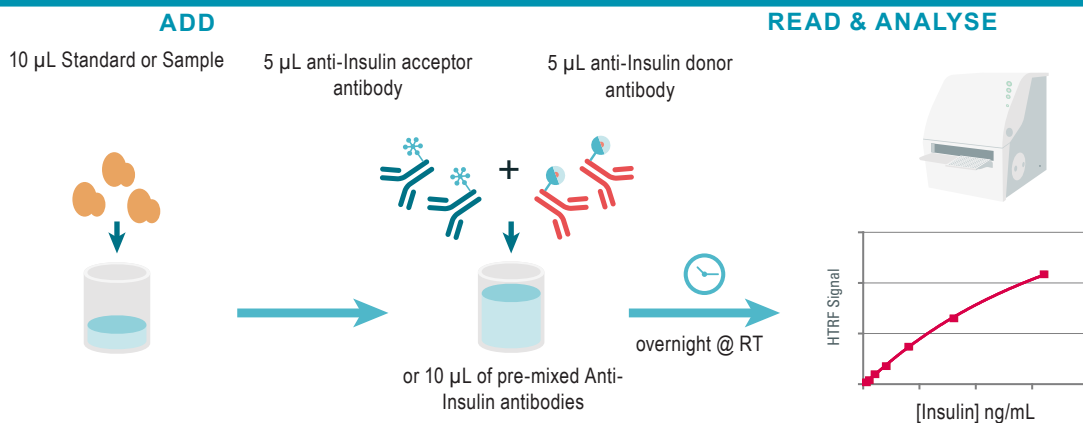


Figure 1: Principle of HTRF Insulin sandwich assay.

MANUAL AT A GLANCE



Make sure to use the set-up for Eu Cryptate.

MATERIALS PROVIDED:

| KIT COMPONENTS | 500 TESTS * CAT # 62IN2PEG | 10,000 TESTS * CAT # 62IN2PEH |
|--------------------------------------|---------------------------------------|---|
| Insulin Standard Lyophilized | 1 vial | 1 vial |
| Insulin Eu Cryptate Antibody | 1 vial Lyophilized | 1 vial Lyophilized |
| Insulin XL665 Antibody | 1 vial Lyophilized | 1 vial Lyophilized |
| Diluent #5 ** 5X | 1 vial 10 mL | 1 vial 100 mL |
| Detection buffer *** ready-to-use | 1 vial 7 mL Detection Buffer #4 | 1 vial 105 mL Detection Buffer #4 |

* When used as advised, the two available kit sizes will provide sufficient reagents for 500 tests and 10,000 tests respectively in 20 µL final volume..

Assay volumes can be adjusted proportionally to run the assay in 96 or 1536 well microplates.

** Medium like cell culture medium can be an alternative to the diluent.

*** The Detection buffer is used to prepare working solutions of acceptor and donor reagents.

PURCHASE SEPARATELY:

- HTRF®-Certified Reader. **Make sure the setup for Eu Cryptate is used.**

For a list of HTRF-compatible readers and set-up recommendations, please visit www.revvity.com

- Small volume (SV) detection microplates - Use white plate only.

For more information about microplate recommendations, please visit our website at: revvity.com

STORAGE AND STABILITY

Store the kit at 2-8°C.

Under proper storage conditions, reagents are stable until the expiry date indicated on the label.

If lyophilized, reconstituted reagents, antibodies, and standard stock solutions may be frozen and thawed only once. To avoid freeze/thaw cycles, it is recommended to dispense remaining stock solutions into disposable plastic vials for storage at -60°C or below .






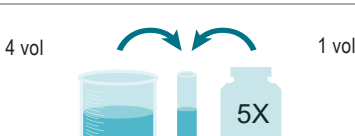
Volume of Insulin Ultra Sensitive standard aliquots should not be under 20 µL.

REAGENT PREPARATION**BEFORE YOU BEGIN:**

- It is very important to prepare reagents in the specified buffers. The use of an incorrect diluent may affect reagent stability and assay results.
- Allow the lyophilized reagents to warm up to room temperature for at least 30 mins before reconstitution
- Before use, allow Diluent and Detection buffer to warm up at room temperature and homogenize them with a vortex.
- It is recommended to filter buffers.
- Antibody solutions must be prepared in individual vials and can be mixed prior to dispensing.
- Insulin standards (for standard curve) must be prepared in diluent or in the same medium as the samples.

TAKE CARE TO PREPARE STOCK AND WORKING SOLUTIONS ACCORDING TO THE DIRECTIONS FOR THE KIT SIZE YOU HAVE PURCHASED.

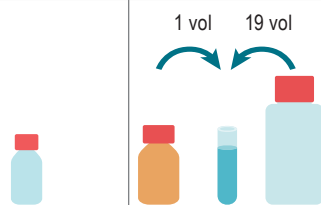
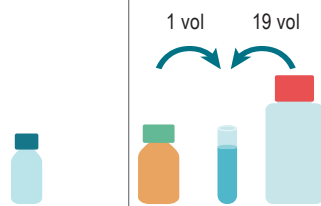
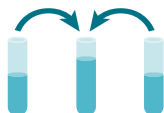
TO PREPARE REAGENT STOCK SOLUTIONS:

| 500 TESTS KIT - 62IN2PEG | | 10,000 TESTS KIT - 62IN2PEH | |
|---|---|---|---|
| Anti-Insulin Eu Cryptate antibody | | | |
| Reconstitute the Insulin Eu Cryptate antibody with 2.5 mL detection buffer #4. Mix gently. This ready-to-use 1X stock solution can be frozen and stored at -60°C or below. |  |  | Reconstitute the Insulin Eu Cryptate antibody with 2.5 mL distilled water. Mix gently. This 20 X stock solution can be frozen and stored at -60°C or below. |
| Anti-Insulin XL665 antibody | | | |
| Reconstitute the Insulin XL665 antibody with 2.5 mL detection buffer #4. Mix gently. This ready-to-use 1X stock solution can be frozen and stored at -60°C or below. |  |  | Reconstitute the Insulin XL665 antibody with 2.5 mL distilled water. Mix gently. This 20 X stock solution can be frozen and stored at -60°C or below. |
| Insulin Standard | | | |
| Reconstitute the Insulin Standard with distilled water in order to obtain a 500 ng/mL stock solution. See instructions on vial label for reconstitution volume. Mix gently after reconstitution. This stock solution can be frozen and stored at -60°C or below. |  | | Reconstitute the Insulin Standard with distilled water in order to obtain a 500 ng/mL stock solution. See instructions on vial label for reconstitution volume. Mix gently after reconstitution. This stock solution can be frozen and stored at -60°C or below. |
| Diluent | | | |
| Dilute 5-fold the 5 X diluent #5 with distilled water: homogenize the 5 X diluent #5 with a vortex and add 1 volume of stock solution in 4 volumes of distilled water (e.g., 1 mL of diluent + 4 mL of distilled water). Mix gently after dilution. This 1X diluent can be frozen and stored at -60°C or below. |  | | Dilute 5-fold the 5 X diluent #5 with distilled water: homogenize the 5 X diluent #5 with a vortex and add 1 volume of stock solution in 4 volumes of distilled water (e.g., 10 mL of diluent + 40 mL of distilled water). Mix gently after dilution. This 1X diluent can be frozen and stored at -60°C or below. |
| Detection buffer | | | |
| The Detection buffer is ready-to-use. | | | The Detection buffer is ready-to-use. |

TO PREPARE ANTIBODY WORKING SOLUTIONS:

Each well requires 5 µL of Insulin-Eu Cryptate Antibody and 5 µL of Insulin-XL665 Antibody.

Prepare the two antibody solutions in separate vials.

| 500 TESTS KIT - 62IN2PEG | 10,000 TESTS KIT - 62IN2PEH | |
|--|---|--|
| Insulin Eu Cryptate antibody | | |
| After reconstitution, the Insulin Eu Cryptate antibody is ready-to-use. |  | Dilute 20-fold the 20 X stock solution (reconstituted reagent) of Insulin Eu Cryptate antibody with the Detection buffer #4: add 1 volume of Eu Cryptate antibody stock solution in 19 volumes of Detection buffer #4 (e.g., 0.5 mL of reconstituted Eu Cryptate antibody stock solution + 9.5 mL of Detection Buffer #4). |
| Insulin XL665 antibody | | |
| After reconstitution, the Insulin XL665-antibody is ready-to-use. |  | Dilute 20-fold the 20 X stock solution (reconstituted reagent) of Insulin XL665 antibody with the Detection buffer #4: add 1 volume of XL665 antibody stock solution in 19 volumes of Detection buffer #4 (e.g., 0.5 mL of reconstituted XL665 antibody stock solution + 9.5 mL of Detection Buffer #4). |
| Antibody mix | | |
| It is possible to pre-mix the two ready-to-use antibody solutions just prior to dispensing the reagents by adding 1 volume of XL665 antibody solution to 1 volume of Cryptate antibody solution (e.g. 1 mL of XL665 antibody + 1 mL of Cryptate antibody). |  | It is possible to pre-mix the two ready-to-use antibody solutions just prior to dispensing the reagents by adding 1 volumes of XL665-antibody solution to 1 volume of Cryptate antibody solution (e.g. 5 mL of XL665-antibody + 5 mL of Cryptate antibody). |

TO PREPARE STANDARD WORKING SOLUTIONS:

- Each well requires 10 μL of standard.
- Dilute the standard stock solution serially with diluent #5 (1X) or with the cell culture medium used to culture the cells (e.g. KRB, KRHB, DMEM, MEM, RPMI+SVF).
- In order to check for a potential interference effect from your own assay buffer when using the assay for the first time, we highly recommend the parallel preparation of a standard curve in your own supplemented cell culture medium and in diluent #5 (1X).
- In order to counteract any standard sticking, we recommend changing tips between each dilution.
- However, if sticking standard is noticed, it is recommended to supplement your assay buffer with 0.05% Tween 20

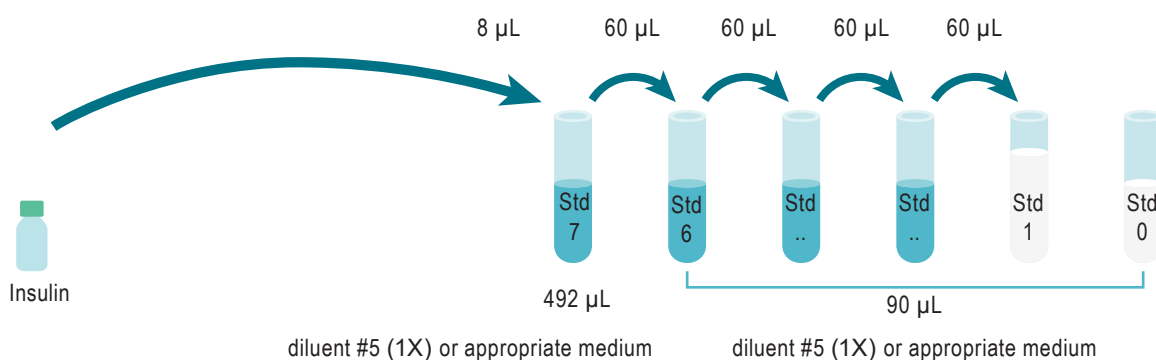
A recommended standard dilution procedure is listed and illustrated below:

Dilute the standard stock solution 62.5-fold with diluent #5 (1X) to prepare high standard (Std 7): e.g. take 8 μL of standard stock solution and add it to 492 μL of diluent #5 (1X). Mix gently.

Use the high standard (Std 7) to prepare the standard curve using 1/2.5 serial dilutions as follows:

- Dispense 90 μL of diluent #5 (1X) in each vial from Std 6 to Std 0.
- Add 60 μL of standard to 90 μL of diluent #5 (1X), mix gently and repeat the 1/2.5 serial dilution to make standard solutions: std6, std5, std4, std3, std2, std1.

This will create 7 standards for the analyte. Std 0 (Negative control) is diluent #5 (1X) or appropriate culture medium alone.








| STANDARD | SERIAL DILUTIONS | INSULIN WORKING SOLUTIONS (ng/mL) |
|-------------------------|--|-----------------------------------|
| Standard Stock solution | Reconstituted lyophilisate | 500 |
| Standard 7 | 8 μL stock solution + 492 μL Diluent #5 (1X) | 8 |
| Standard 6 | 60 μL standard 7 + 90 μL Diluent #5 (1X) | 3.2 |
| Standard 5 | 60 μL standard 6 + 90 μL Diluent #5 (1X) | 1.28 |
| Standard 4 | 60 μL standard 5 + 90 μL Diluent #5 (1X) | 0.5 |
| Standard 3 | 60 μL standard 4 + 90 μL Diluent #5 (1X) | 0.2 |
| Standard 2 | 60 μL standard 3 + 90 μL Diluent #5 (1X) | 0.08 |
| Standard 1 | 60 μL standard 2 + 90 μL Diluent #5 (1X) | 0.03 |
| Standard 0 | 90 μL Diluent #5 (1X) | 0 |

TO PREPARE SAMPLES:

- Each well requires 10 μ L of sample.
- Just after their collection, put the samples at 4°C and test them immediately. For later use, samples should be dispensed into disposable plastic vials and stored at -60°C or below. Avoid multiple freeze/thaw cycles.
- In order to measure the total Insulin content in cell or islet lysates, we recommend an acid-ethanol extraction step
- Samples with a concentration above the highest standard (Std 7) must be diluted diluent #5 (1X)
- Before measurement, cell or islet lysates must be diluted at least 100-fold in the used secretion assay buffer

ASSAY MANUAL

| | | Standard (Std 0 - Std 7) | Samples |
|--------|---|--|--|
| Step 1 |  | Dispense 10 μ L of each Insulin standard (Std 0 - Std 7) into each standard well | Dispense 10 μ L of each sample into each sample well |
| Step 2 |  | Add 5 μ L of Insulin XL665 antibody working solution to all wells | |
| Step 3 |  | Add 5 μ L of Insulin Eu Cryptate antibody working solution to all wells | |
| Step 4 |  | Seal the plate and incubate overnight @ RT | |
| Step 5 |  | Remove the plate sealer and read on an HTRF® compatible reader | |

| | 1 | 2 | 3 | 4 | 5 | 6 |
|---|--|----------------|----------------|--|----------------|----------------|
| A | 10 µL Std 0 (Negative control) 5 µL Insulin-XL665 5 µL Insulin-Eu Cryptate | Repeat Well A1 | Repeat Well A1 | 10 µL Sample 1 5 µL Insulin-XL665 5 µL Insulin-Eu Cryptate | Repeat Well A4 | Repeat Well A4 |
| B | 10 µL Std 1 5 µL Insulin-XL665 5 µL Insulin-Eu Cryptate | Repeat Well B1 | Repeat Well B1 | 10 µL Sample 2 5 µL Insulin-XL665 5 µL Insulin-Eu Cryptate | Repeat Well B4 | Repeat Well B4 |
| C | 10 µL Std 2 5 µL Insulin-XL665 5 µL Insulin-Eu Cryptate | Repeat Well C1 | Repeat Well C1 | 10 µL Sample 3 5 µL Insulin-XL665 5 µL Insulin-Eu Cryptate | Repeat Well C4 | Repeat Well C4 |
| D | 10 µL Std ... 5 µL Insulin-XL665 5 µL Insulin-Eu Cryptate | Repeat Well D1 | Repeat Well D1 | 10 µL Sample ... 5 µL Insulin-XL665 5 µL Insulin-Eu Cryptate | Repeat Well D4 | Repeat Well D4 |
| E | 10 µL Std ... 5 µL Insulin-XL665 5 µL Insulin-Eu Cryptate | Repeat Well E1 | Repeat Well E1 | 10 µL Sample ... 5 µL Insulin-XL665 5 µL Insulin-Eu Cryptate | Repeat Well E4 | Repeat Well E4 |
| F | 10 µL Std ... 5 µL Insulin-XL665 5 µL Insulin-Eu Cryptate | Repeat Well F1 | Repeat Well F1 | 10 µL Sample ... 5 µL Insulin-XL665 5 µL Insulin-Eu Cryptate | Repeat Well F4 | Repeat Well F4 |
| G | 10 µL Std ... 5 µL Insulin-XL665 5 µL Insulin-Eu Cryptate | Repeat Well G1 | Repeat Well G1 | 10 µL Sample ... 5 µL Insulin-XL665 5 µL Insulin-Eu Cryptate | Repeat Well G4 | Repeat Well G4 |
| H | 10 µL Std ... 5 µL Insulin-XL665 5 µL Insulin-Eu Cryptate | Repeat Well H1 | Repeat Well H1 | 10 µL Sample ... 5 µL Insulin-XL665 5 µL Insulin-Eu Cryptate | Repeat Well H4 | Repeat Well H4 |

DATA REDUCTION

1. Calculate the ratio of the acceptor and donor emission signals for each individual well.

$$\text{Ratio} = \frac{\text{Signal 665 nm}}{\text{Signal 620 nm}} \times 10^4$$

2. Calculate the % CVs. The mean and standard deviation can then be worked out from ratio replicates.

$$\text{CV (\%)} = \frac{\text{Standard deviation}}{\text{Mean Ratio}} \times 100$$

3. Calculate the % delta F which reflects the signal to background of the assay. The negative control (Standard 0) plays the role of an internal assay control. Delta F is used for the comparison of day to day runs of the same assay.

$$\text{delta F (\%)} = \frac{\text{Ratio Standard or sample} - \text{Ratio Negative Control}}{\text{Ratio Negative Control}} \times 100$$

For more information about data reduction, please visit www.revvity.com

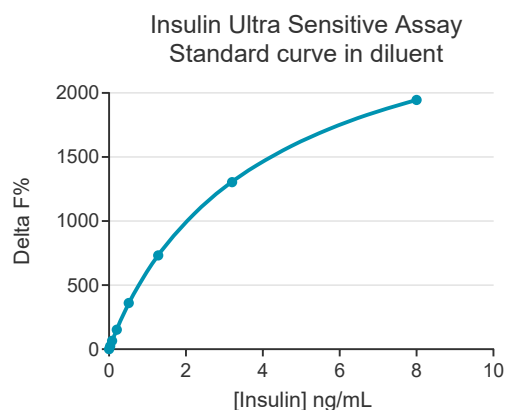
RESULTS

This data must not be substituted for the data obtained in the laboratory and should be considered only as an example.

Results may vary from one HTRF® compatible reader to another.

Standard curve fitting with the 4 Parameter Logistic (4PL) model:

| | Ratio ⁽¹⁾ | CV ⁽²⁾ | Delta F% ⁽³⁾ |
|-------------------------------|----------------------|-------------------|-------------------------|
| Standard 0 - Negative control | 626 | 0.2% | - |
| Standard 1 - 0.033 ng/mL | 779 | 0.9% | 24% |
| Standard 2 - 0.082 ng/mL | 1,035 | 2% | 65% |
| Standard 3 - 0.205 ng/mL | 1,628 | 2.3% | 160% |
| Standard 4 - 0.512 ng/mL | 2,914 | 5.2% | 365% |
| Standard 5 - 1.28 ng/mL | 5,741 | 1.5% | 817% |
| Standard 6 - 3.2 ng/mL | 11,335 | 2.4% | 1,710% |
| Standard 7 - 8 ng/mL | 17,472 | 3.4% | 2,690% |



ANALYTICAL CHARACTERISTICS

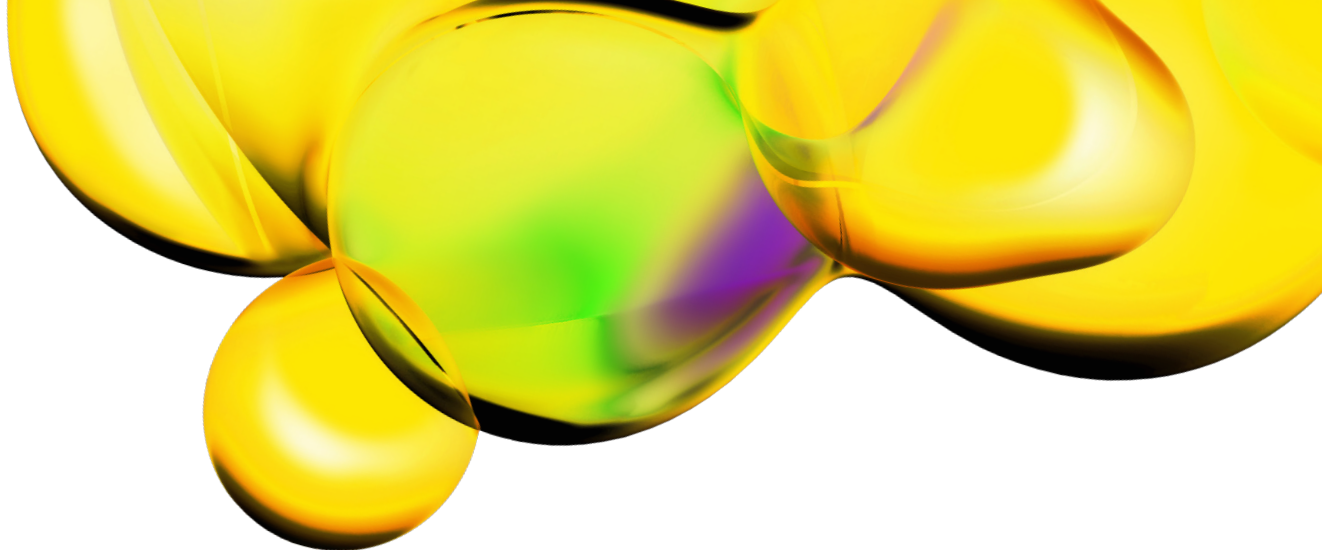
ASSAY PERFORMANCES AND CALIBRATION

| | |
|---|--|
| Calibration | NIBSC International standard 66/304 - 1 ng Insulin HTRF \longleftrightarrow 1 ng Insulin NIBSC 66/304 (i.e.: 2.3×10^{-5} IU) |
| Assay range | 0.033 to 8 ng/mL |
| Limit of detection (LoD*) = Std 0 Mean + 2 SD | 0.005 ng/mL |
| incubation time | Overnight at RT |

*The LoD was calculated from data obtained with the PHERAstar Plus reader (flash lamp excitation) after overnight incubation. This may vary from one HTRFcompatible reader to another.

This product contains material of biologic origin. Use for research purposes only. Do not use in humans or for diagnostic purposes. The purchaser assumes all risk and responsibility concerning reception, handling and storage.

The use of the cell line will be done with appropriate safety and handling precautions to minimize health and environmental impact.



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF

Manufactured by Cisbio Bioassays - 172 rue de la Cèze - Parc Marcel Boiteux - 30200 Codolet - FRANCE

www.revvity.com

revvity

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2025, Revvity, Inc. All rights reserved.