



## LANCE *Ultra* c-Myc (Human) Detection Kit

**Product number:** TRF1111 C/M

Research Use Only. Not for use in diagnostic procedures.

### Product Information

**Application:** This kit is designed for the quantitative determination of c-Myc in media using a homogeneous LANCE *Ultra* assay (no wash steps).

**Sensitivity:** Lower Detection Limit (LDL): 46.1 nM  
Lower Limit of Quantification (LLOQ): 235.8 nM  
EC<sub>50</sub>: 1 346.4 nM

**Dynamic range:** Kit designed to detect c-Myc between: 46.1– 300,000 nM (Figure 1).

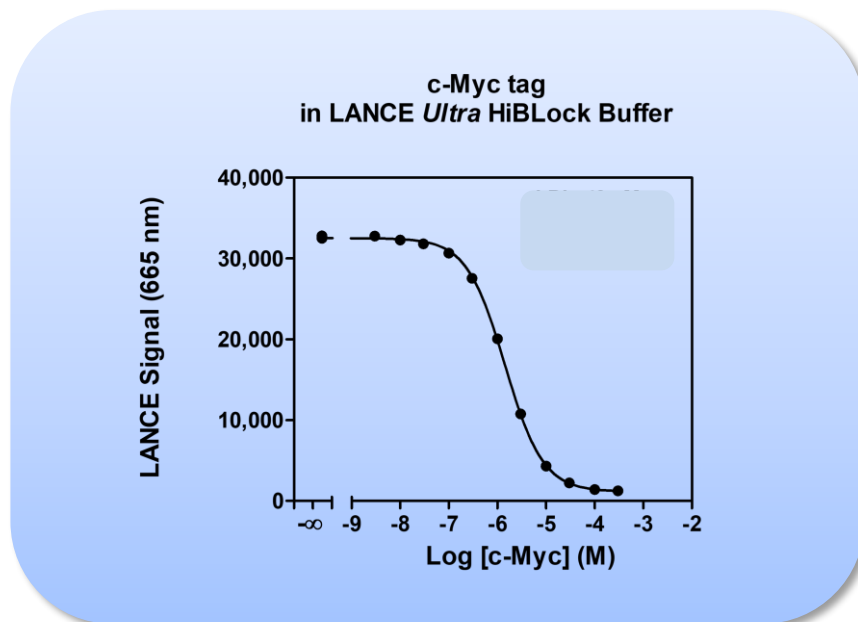


Figure 1. Typical sensitivity curves in *Ultra* HiBlock Buffer. The data was generated using a white OptiPlate™ -384 microplate and the EnVision Plate Reader equipped with TR-FRET option and a TRF laser. Performance on other plate readers may vary.

**Storage:** Store kit in the dark at +4°C. Store analyte at -20°C upon receipt. Store reconstituted analyte at - 20°C.

**Stability:** This kit is stable for at least 12 months from the manufacturing date when stored in its original packaging and the recommended storage conditions.

## Analyte of Interest

The c-myc polypeptide tag is derived from the c-myc gene. The tag can be used during affinity chromatography and to separate wild-type protein from recombinant protein. The c-myc tag can be used in indirect detection assays for which there is no antibody for the target.

## Description of the LANCE *Ultra* Assay

LANCE® and LANCE® (Lanthanide chelate excite) *Ultra* are our TR-FRET (time-resolved fluorescence resonance energy transfer), homogeneous (no wash) technologies. One antibody of interest is labeled with an acceptor fluorophore [*ULight*™ dye]. A “tracer” analyte is biotinylated and binds Streptavidin labeled with donor Eu-chelate. Upon excitation at 320 or 340 nm, energy can be transferred from the donor Europium chelate to the acceptor fluorophore if sufficiently close for FRET (~10 nm). This results in the emission of light at 665 nm. In the presence of sufficient non-biotinylated analyte this emission will be diminished.

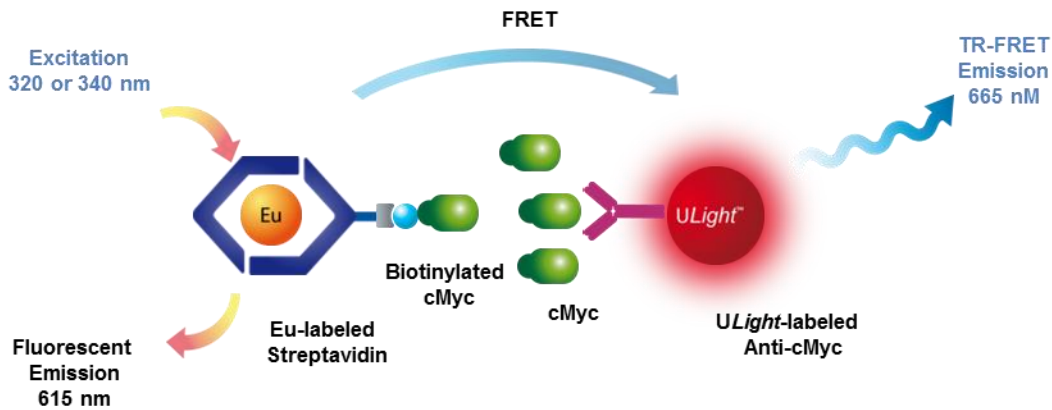


Figure 2. LANCE competition assay principle.

## Precautions

- All blood components and biological materials should be handled as potentially hazardous.
- Some analytes are present in saliva. Take precautionary measures to avoid contamination of the reagent solutions.

## Kit Content: Reagents and Materials

Kit components	TRF1111C (500 assay points <sup>***</sup> )	TRF1111M (10 000 assay points <sup>***</sup> )
LANCE <i>Ultra ULight</i> -labeled Anti- c-Myc Antibody stored in TSA buffer, 0.1% BSA	60 µL @ 500 nM (1 brown tube, blue cap)	1200 µL @ 500 nM (1 brown tube, green cap)
Biotinylated c-Myc-Tracer	25 µL @ 100 µM (1 tube, black cap)	25 µL @ 10 µM (1 tube, black cap)
c-Myc Analyte* standard stored in 1% Acetic Acid	100 µL @ 10 mM (1 tube, <u>clear</u> cap)	100 µL @ 10 mM (1 tube, <u>clear</u> cap)
LANCE Eu-W1024-labeled Streptavidin	20 µL @ 8.3 µM (1 clear tube, yellow cap)	20 µL @ 1.3 µM (1 clear tube, orange cap)
<i>Ultra</i> HiBlock Buffer (5X)**	2 mL, 1 small bottle	100 mL, 1 large bottle

\* One vial contains an amount of c-Myc sufficient for performing 10 standard curves. Aliquot and store unused analyte at -20 °C. Additional vials can be ordered separately (cat # TRF1111S).

\*\* Extra buffer can be ordered separately (cat # TRF1011C: 10 mL, cat # TRF1011F: 100 mL). 5X *Ultra* HiBlock Buffer may appear cloudy, especially after storage at cold temperature. Agitate and/or stir at room temperature to redissolve prior to dilution.

\*\*\* The number of assay points is based on an assay volume of 20 µL in 384-well assay plates using the kit components at the recommended concentrations.

Sodium azide should **not** be added to the stock reagents. High concentrations of sodium azide (> 0.001 % final in the assay) might decrease the signal.

### Specific additional required reagents and materials:

The following materials are recommended:

Item	Suggested source	Catalog #
TopSeal-A PLUS Adhesive Sealing Film	Revvity Inc.	6050185
EnVision Multilabel Plate Reader equipped with TR-FRET option and a TRF laser	Revvity Inc.	-

## Recommendations

### General recommendations:

- The volume indicated on each tube is guaranteed for single pipetting. Multiple pipetting of the reagents may reduce the theoretical amount left in the tube.
- Centrifuge all tubes (including analyte) before use to improve recovery of content (2000g, 10-15 sec).
- Re-suspend all reagents by vortexing before use.
- Use Milli-Q® grade H<sub>2</sub>O (18 MΩ•cm) to dilute Buffer.
- When diluting the standard or samples, change tips between each standard or sample dilution. When loading reagents in the assay microplate, change tips between each standard or sample addition and after each set of reagents.
- When reagents are added to the microplate, make sure the liquids are at the bottom of the well.
- Small volumes may be prone to evaporation. It is recommended to cover microplates with TopSeal-A Adhesive Sealing Films to reduce evaporation during incubation. LANCE Ultra TR-FRET assays cannot be read with the TopSeal-A Film attached. Please remove before reading.
- LANCE signal is detected using a VICTOR X, ViewLux, EnVision or EnSpire Multilabel Reader equipped with the TR-FRET. Use an excitation wavelength of 320 or 340 nm to excite the LANCE Europium chelate. We recommend you read this assay in dual emission mode, detecting both the emission from the Europium donor fluorophore at 615 nm, and the acceptor fluorophore (at 665 nm for *ULight* dye). The raw FRET signal at 665 nm can be used to process your data.
- Signal will vary with temperature and incubation time. For consistent results, identical incubation times and temperature should be used for each plate.
- The standard curves shown in this technical data sheet are provided for information only. A standard curve must be generated for each experiment. The standard curve should be performed in *Ultra* HiBlock Buffer

## Assay Procedure

IMPORTANT: PLEASE READ THE RECOMMENDATIONS BELOW BEFORE USE

- The manual described below is an **example** for generating one standard curve in a 20  $\mu\text{L}$  final assay volume (48 wells, triplicate determinations) and 452 samples. The manuals also include testing samples in 384 well plates. If different amounts of samples are tested, the volumes of all reagents must be adjusted accordingly, as shown in the table below. \*\*\* These calculations do not include excess reagents to account for losses during transfer of solutions or dead volumes.
- The standard dilution manual is provided for information only. As needed, the number of replicates or the range of concentrations covered can be modified.
- Use of four background points in triplicate (12 wells) is recommended when LDL/LLOQ is calculated. One background point in triplicate (3 wells) can be used when LDL/LLOQ is not calculated.

Format	# of data points	Volume			Plate recommendation
		Final	Sample	Eu-Antibody/ULight-Antibody MIX	
TRF1111C	250	40 $\mu\text{L}$	30 $\mu\text{L}$	10 $\mu\text{L}$	White OptiPlate-96 (cat # 6005290) White $\frac{1}{2}$ AreaPlate-96 (cat # 6005560)
	500	20 $\mu\text{L}$	15 $\mu\text{L}$	5 $\mu\text{L}$	White $\frac{1}{2}$ AreaPlate-96 (cat # 6005560) White OptiPlate-384 (cat # 6007290)
	1 250	8 $\mu\text{L}$	6 $\mu\text{L}$	2 $\mu\text{L}$	ProxiPlate™-384 Plus (cat # 6008280) White OptiPlate-384 (cat # 6007290)
	2 500	4 $\mu\text{L}$	3 $\mu\text{L}$	1 $\mu\text{L}$	White OptiPlate-1536 (cat # 6004290)
TRF1111M	10 000	20 $\mu\text{L}$	15 $\mu\text{L}$	5 $\mu\text{L}$	White $\frac{1}{2}$ AreaPlate-96 (cat # 6005560) White OptiPlate-384 (cat # 6007290)
	25 000	8 $\mu\text{L}$	6 $\mu\text{L}$	2 $\mu\text{L}$	ProxiPlate-384 Plus (cat # 6008280) White OptiPlate-384 (cat # 6007290)
	50 000	4 $\mu\text{L}$	3 $\mu\text{L}$	1 $\mu\text{L}$	White OptiPlate-1536 (cat # 6004290)

**General Manual (1-step manual):** Dilute standards, samples, and assay components in 1X *Ultra* HiBlock Buffer.

**Each manual described below is designed for 500 assay points including one standard curve (48 wells) and samples (452 wells).**

**Standard Preparation:**1) Preparation of 1X *Ultra* HiBlock Buffer:

- a. Add 2 mL of 5X *Ultra* HiBlock Buffer to 8 mL H<sub>2</sub>O.

2) Preparation of c-Myc analyte standard dilutions:

- a. c-Myc analyte is provided at 10 mM in 1% Acetic Acid. Prepare standard dilutions as follows (change tip between each standard dilution):
- b. **Tube A may become cloudy after addition of 10 mM stock. Exclude point from analysis.**

Tube	Vol. of c-Myc (µL)	Vol. of diluent (µL) *	[c-Myc] in standard curve
			(M in 5 µL)
A	10 µL of 10 mM stock c-Myc	90	1.0E-3
B	30 µL of tube A	70	3.0E-4
C	30 µL of tube B	60	1.0E-4
D	30 µL of tube C	70	3.0E-5
E	30 µL of tube D	60	1.0E-5
F	30 µL of tube E	70	3.0E-6
G	30 µL of tube F	60	1.0E-6
H	30 µL of tube G	70	3.0E-7
I	30 µL of tube H	60	1.0E-7
J	30 µL of tube I	70	3.0E-8
K	30 µL of tube J	60	1.0E-8
L	30 µL of tube K	70	3.0E-9
M ** (background)	0	100	0
N ** (background)	0	100	0
O ** (background)	0	100	0
P ** (background)	0	100	0

\* At low concentrations of analyte, a significant amount of analyte can bind to the vial. Therefore, load the analyte standard dilutions in the assay microplate within 60 minutes of preparation.

\*\* Four background points in triplicate (12 wells) are used when LDL is calculated. If LDL does not need to be calculated, one background point in triplicate can be used (3 wells).

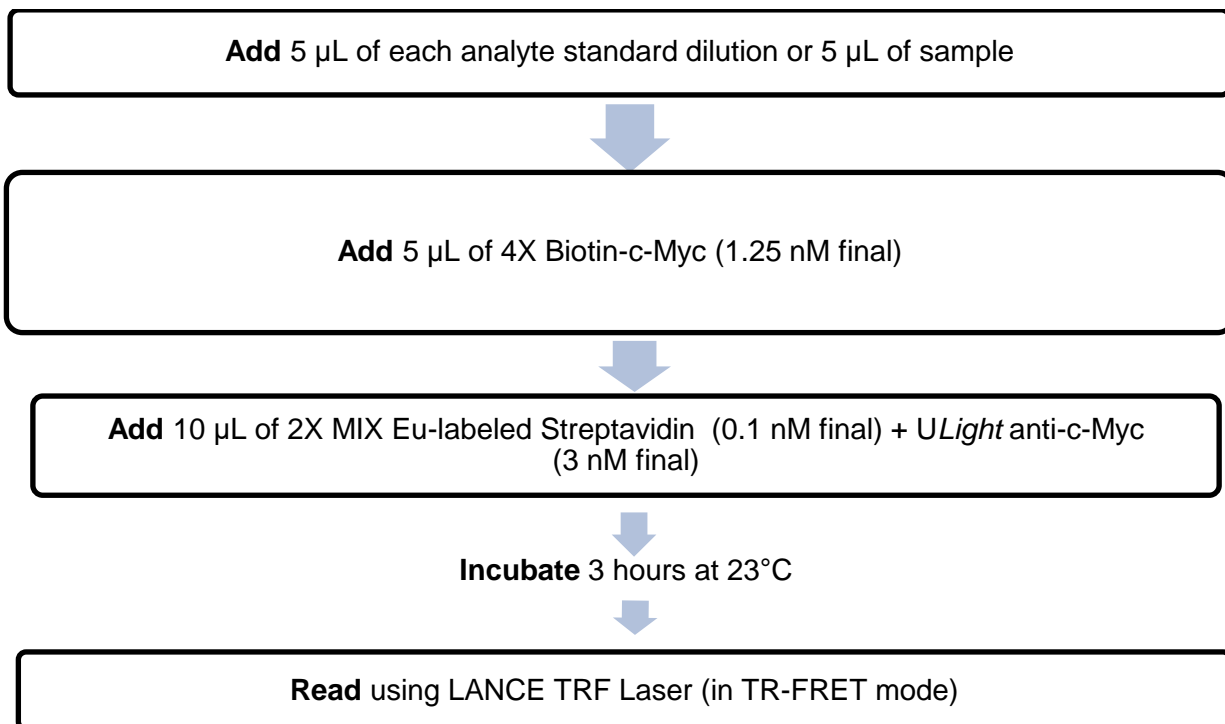
3) Preparation of 4X biotin-c-Myc (tracer) (5 nM):

- a. Prepare a 500X biotin-cMyc Tracer:
  - i. Add 3.125 µL of 100 µM stock to 496.875 µL *Ultra* HiBlock Buffer
- b. Add 4 µL of 500 biotin-c-Myc to 496 µL of *Ultra* HiBlock Buffer.
- c. Prepare just before use.

4) Preparation of 2X MIX Eu-labeled Streptavidin (0.2 nM) + *ULight* labeled anti-c-Myc Antibody (6 nM):

- a. Prepare 500X Eu-SA:
  - i. Add 3 µL of 8.3 µM Eu-SA to 497 µL *Ultra* HiBlock Buffer
- b. Add 12 µL of 500 nM *ULight* labeled anti-c-Myc Antibody and 4 µL of 500X Europium labeled Streptavidin to 984 µL of *Ultra* HiBlock Buffer.
- c. Prepare just before use.

5) In a white OptiPlate (384 wells):



**Important:** LANCE signal is detected using an EnVision Multilabel Reader equipped with the TR-FRET. Use an excitation wavelength of 320 or 340 nm to excite the LANCE Europium chelate. We recommend you read this assay in dual emission mode, detecting both the emission from the Europium donor fluorophore at 615 nm, and the acceptor fluorophore (at 665 nm for *ULight* dye).

## Data Analysis

- Calculate the average count value for the background wells.
- Generate a standard curve by plotting the LANCE counts versus the concentration of analyte. A log scale can be used for either or both axes. No additional data transformation is required.
- Analyze data according to a nonlinear regression using the 4-parameter logistic equation (sigmoidal dose-response curve with variable slope) and a  $1/Y^2$  data weighting (the values at maximal concentrations of analyte after the hook point should be removed for correct analysis).
- The LDL is calculated by interpolating the average background counts (12 wells without analyte) + 2 x standard deviation value (average background counts + (2xSD)) on the standard curve.
- The LLOQ as measured here is calculated by interpolating the average background counts (12 wells without analyte) + 10 x standard deviation value (average background counts + (10xSD)) on the standard curve. Alternatively, the true LLOQ can be determined by spiking known concentrations of analyte in the matrix and measuring the percent recovery, and then determining the minimal amount of spiked analyte that can be quantified within a given limit (usually +/- 20% or 30% of the real concentration).
- Read from the standard curve the concentration of analyte contained in the samples.
- If samples have been diluted, the concentration read from the standard curve must be multiplied by the dilution factor.

## Assay Performance Characteristics

LANCE Ultra assay performance described below was determined using the 1 step manual.

### Assay Sensitivity

The LDL and LLOQ were calculated as described above. The values correspond to the lowest concentration of analyte that can be detected in a volume of 15  $\mu$ L using the recommended assay conditions.

LDL (nM)	LLOQ (nM)	Buffer	# of experiments
46	238	<i>Ultra</i> HiBlock	9
37	251	DMEM + 0.1% BSA	6

### Assay Precision:

The following assay precision data were calculated from the three independent assays using two different kit lots. In each lot, the analytes were prepared in *Ultra* HiBlock Buffer. Each assay consisted of one standard curve comprising 12 data points in triplicate and 12 background wells containing no analyte. The assays were performed in a 384-well format using *Ultra* HiBlock Buffer.

#### Intra-assay precision:

The intra-assay precision was determined using 3 independent experiments for a total of 16 independent determinations in triplicate. CV% were calculated for each individual experiment then averaged. Shown is the average intra-experimental CV%.

c-Myc (CV%)	Buffer
3	<i>Ultra</i> HiBlock
3	DMEM + 0.1% BSA

#### Inter-assay precision:

The inter-assay precision was determined using the data across 3 independent experiments with 16 measurements in triplicate. CV% was calculated by comparing the same measurement in each experiment. The CV% for all 16 measurements were then averaged. Shown is the inter-experimental CV%.

c-Myc (CV%)	Buffer
11	<i>Ultra</i> HiBlock
4	DMEM +0.1% BSA



### **Spike Recovery:**

In four experiments, three known concentrations of c-Myc were spiked into 2 separate media and performed triplicate. The spiked samples were referenced to the c-Myc analyte curve produced in the corresponding media.

Spiked c-Myc ( $\mu\text{M}$ )	% Recovery	
	<i>Ultra HiBlock Buffer</i>	DMEM + 0.1% BSA
10	98	87
3	96	96
1	96	90

### **Troubleshooting Guide**

You will find detailed recommendations for common situations you might encounter with your LANCE *Ultra* Assay kit at:

[www.revivity.com](http://www.revivity.com)

**FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.**



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

[www.revvity.com](http://www.revvity.com)

revvity

**Revvity, Inc.**  
940 Winter Street  
Waltham, MA 02451 USA  
[www.revvity.com](http://www.revvity.com)

For a complete listing of our global offices, visit [www.revvity.com](http://www.revvity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.