

LANCE *Ultra* EGF and EGFR (Human) Binding Kit

Product number: TRF1366 C/M

Caution: For Laboratory Use. A research product for research purposes only.

Product Information

Application: This kit is designed for the detection of binding activity between Human EGF and EGFR using a homogeneous LANCE *Ultra* assay (no wash steps). This assay can facilitate the design and development of therapeutics which competitively inhibit EGF/EGFR interaction.

Sensitivity: $K_{d(app)}$: 0.25 nM (average) using 3 nM EGFR

Signal to background ratio: 10 (average) using 3 nM EGFR

Storage: The kit components must be stored at +4 °C in the dark.

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.

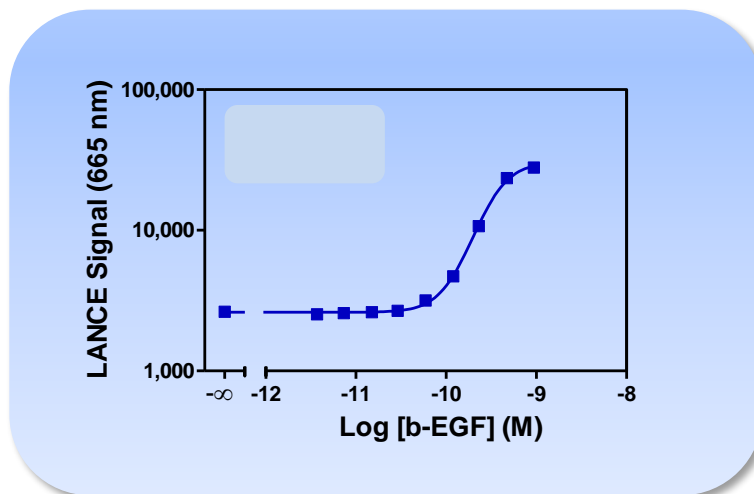


Figure 1. Typical sensitivity curves in *Ultra* HiBlock Buffer. The data was generated using a white OptiPlate™ -384 microplate and the EnVision Plate Reader equipped with TR-FRET option and a TRF laser. Performance on other plate readers may vary.

Analyte of Interest

Epidermal Growth Factor Receptor (EGFR, Her1, ErbB1) is a 134 kDa cell surface receptor, part of a four member subfamily of receptor tyrosine kinases (erbB1, erbB2, erbB3, and erbB4). EGFR binds to a family of proteins called the epidermal growth factors (primarily EGF), which, upon association induces dimerization of the receptor and initiates signal transduction for the promotion of cell survival and growth. Mutations that cause aberrant overexpression of EGF and/or EGFR have been linked with many different cancers such as non-small-cell lung and colon cancers. Drugs are being developed as a means to better treat these types cancer.

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. Streptavidin is labeled with a donor fluorophore (a LANCE Europium chelate) which binds to biotinylated EGF and the EGFR-Fc is recognized by anti-human IgG labeled with an acceptor fluorophore [ULight™ dye]. 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

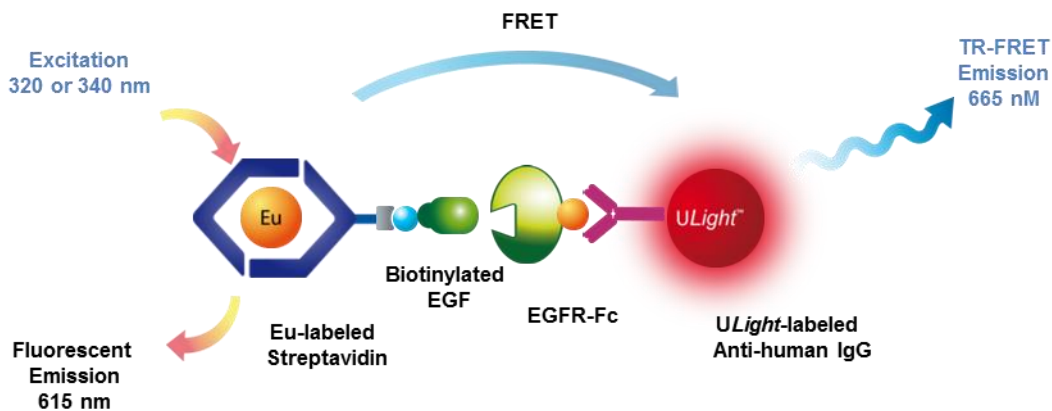


Figure 2. LANCE *Ultra* Assay Principle.

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	TRF1366C (500 assay points)	TRF1366M (10 000 assay points)
LANCE Eu-W1024 labeled Streptavidin stored in TSA buffer, 0.1%BSA	18 µL @ 10 µg/mL (1 clear tube, yellow cap)	360 µL @ 10 µg/mL (1 clear tube, orange cap)
LANCE <i>Ultra ULight</i> -labeled Anti-hIgG stored in TSA buffer, 0.1% BSA	200 µL @ 500 nM (1 brown tube, blue cap)	4 X 1 mL @ 500 nM (4 brown tubes, green caps)
Human EGF (Biotinylated) lyophilized	0.3 µg (1 tube, <u>clear</u> cap)	0.3 µg (8 tubes, <u>clear</u> cap)
Human EGFR -Fc lyophilized	9.56 µg (1 tube, <u>clear</u> cap)	9.56 µg (6 tubes, <u>clear</u> cap)
<i>Ultra</i> HiBlock Buffer (5X)	2 mL, 1 small bottle	100 mL, 1 large bottle

* Reconstitute biotinylated EGF and EGFR in 100 and 200 µL Milli-Q® grade H₂O respectively. The reconstituted proteins should be used within 60 minutes or aliquoted into screw-capped polypropylene vials and stored at -20°C for further experiments. Avoid multiple freeze-thaw cycles.

** 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 Optiplates or 20 µL in 384 well ProxiPlates 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 LANCE *Ultra* 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

- 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 lyophilized analyte) before use to improve recovery of content (2000g, 10-15 sec).
- Re-suspend all reagents by vortexing before use.
- Use Milli-Q® grade H2O (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

Competition Assay Procedure

Format	# of data points	Final	Volume				Plate recommendation
			Inhibitor Or Antibody	EGFR	Biotinylated EGF	ULight anti-IgG + Eu-SA	
TRF1366C	250	40 µL	10 µL	10 µL	10 µL	10 µL	White OptiPlate-96 (cat # 6005290) White ½ AreaPlate-96 (cat # 6005560)
	500	20 µL	5 µL	5 µL	5 µL	5 µL	White ½ AreaPlate-96 (cat # 6005560) White OptiPlate-384 (cat # 6007290)
	1 250	8 µL	2 µL	2 µL	2 µL	2 µL	ProxiPlate™-384 Plus (cat # 6008280) White OptiPlate-384 (cat # 6007290)
	2 500	4 µL	1 µL	1 µL	1 µL	1 µL	White OptiPlate-1536 (cat # 6004290)

TRF1366M	10000	20 µL	10 µL	10 µL	10 µL	10 µL	White ½ AreaPlate-96 (cat # 6005560) White OptiPlate-384 (cat # 6007290)
	25 000	8 µL	2 µL	2 µL	2 µL	2 µL	ProxiPlate-384 Plus (cat # 6008280) White OptiPlate-384 (cat # 6007290)
	50 000	4 µL	1 µL	1 µL	1 µL	1 µL	White OptiPlate-1536 (cat # 6004290)

Example: Competition using target and non-target antibodies

- The manual described below is an **example** for generating 500 wells, triplicate determinations, including three inhibition curves in a 20 µL final assay volume. If a different amount of samples are tested, the volumes of all reagents have to be adjusted accordingly, as shown in the table below. These calculations do not include excess reagent to account for losses during transfer of solutions or dead volumes.
- The dilution manual is provided for information only. As needed, the number of replicates or the range of concentrations covered can be modified.

The following reagents were used for the application example shown below:

Item (Ab)	Supplier	Catalog number
Anti-human EGFR antibody	R & D Systems	AF231

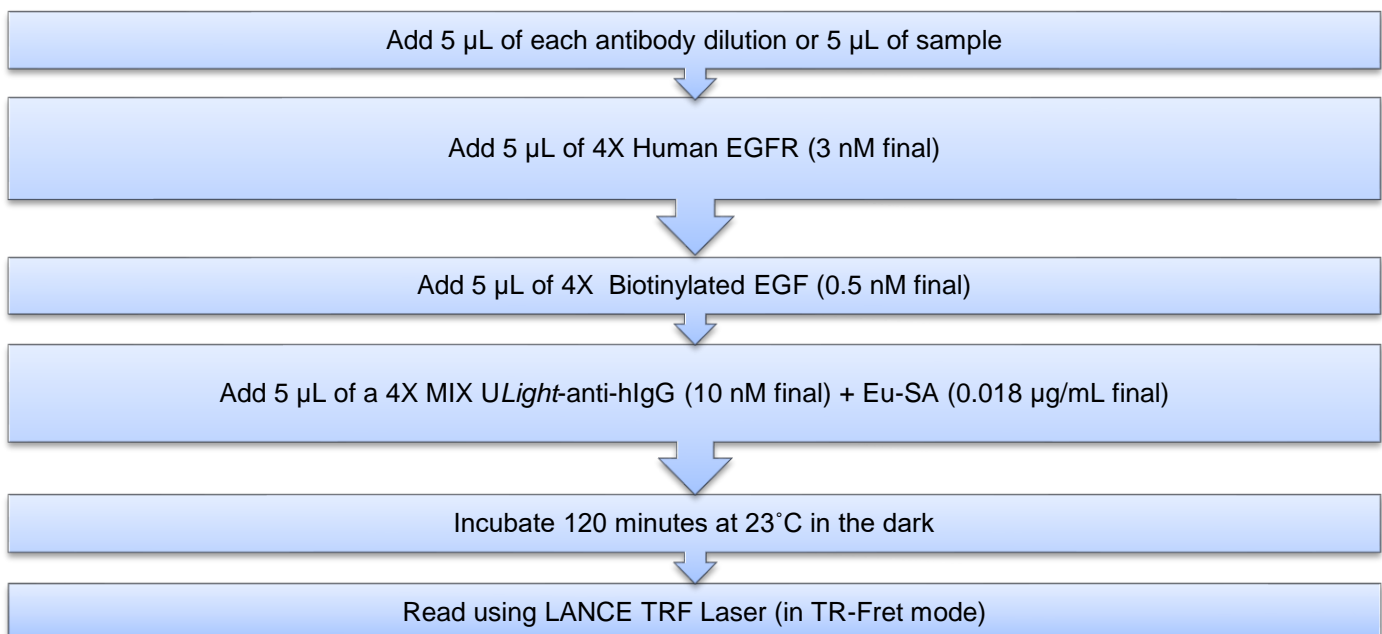
1 step manual (1 Incubation steps) described as below:

- Preparation of 1X *Ultra* HiBlock Buffer:
Add 2 mL of 5X *Ultra* HiBlock Buffer to 8 mL H₂O.
- Preparation of serial dilutions of 4X anti-human EGFR antibody in 1X *Ultra* HiBlock Buffer as follows:

Tube	Volume of Antibody	Volume of 1X buffer	[Ab] (g/mL) (1X) final concentration
A	Add 30 µL of 1E-3 g/mL stock	45	1.00E-04
B	30 µL of tube A	70	3.00E-05
C	30 µL of tube B	60	1.00E-05
D	30 µL of tube C	70	3.00E-06
E	30 µL of tube D	60	1.00E-06
F	30 µL of tube E	70	3.00E-07
G	30 µL of tube F	60	1.00E-07
H	30 µL of tube G	70	3.00E-08

I	30 μ L of tube H	60	1.00E-08
J	30 μ L of tube I	70	3.00E-09
K	30 μ L of tube J	60	1.00E-09
L	30 μ L of tube K	70	3.00E-10
M-Max Signal	0	100 μ L	0
N-Max Signal	0	100 μ L	0
O-Max Signal	0	100 μ L	0
P-Max Signal	0	100 μ L	0

- 3) Preparation of 4X Mix of *ULight* labeled Anti-Human IgG Fc-specific Antibody (40 nM) and Eu-W1024 labeled-Streptavidin (0.072 μ g/mL):
- Add 18 μ L of Eu-W1024 Streptavidin at 10 μ g/mL and 200 μ L of *ULight* anti Human IgG at 500 nM to 2282 μ L of 1X *Ultra* HiBlock Buffer
 - Prepare just before use.
- 4) Preparation of 4X biotinylated EGF(2 nM):
- Prepare just before use.
 - Reconstitute lyophilized biotinylated EGF (0.3 μ g) in 100 μ L H₂O to make 500 nM
 - Add 10 μ L of 500 nM biotinylated EGF to 2490 μ L 1X *Ultra* HiBlock Buffer.
- 5) Preparation of 4X EGFR Fc-tagged (12 nM):
- Reconstitute lyophilized EGFR Fc-tagged (9.56 μ g) in 200 μ L H₂O to make 500 nM stock concentration.
 - Add 60 μ L of 500 nM EGFR Fc-tagged to 2440 μ L of 1X *Ultra* HiBlock Buffer.
 - Prepare just before use.
- 6) 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).

Sample Results:

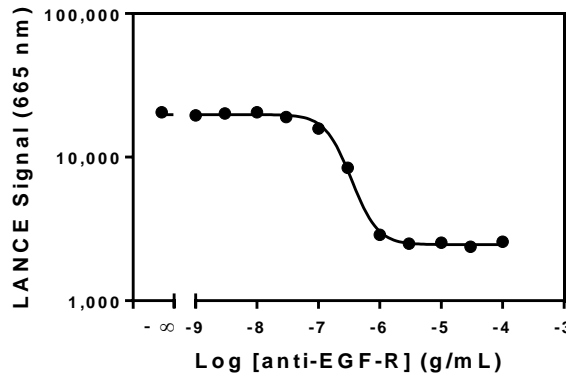


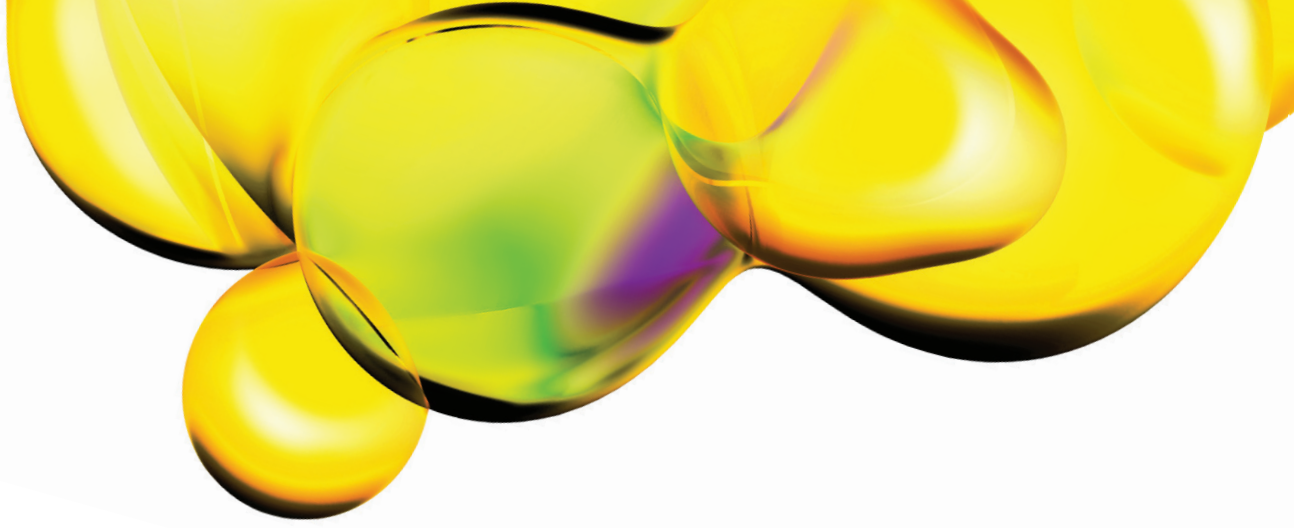
Figure 3. Antibody Inhibition curve using EGFR antibody. Anti-Human EGFR (calculated IC₅₀ of 215 nM)

Troubleshooting Guide

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

www.revvy.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

revvity

Revvity, Inc.
940 Winter Street
Waltham, MA 02451 USA
www.revvity.com

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2023, Revvity, Inc. All rights reserved.