

HTRF® Research Reagents

HTRF EGFR total kit - 10K pts

Product Number : 64NG1PEH

Lot Number : 12G

Expiration date : 2027-01-20

Document version :

Reagents :	Lot n°
HTRF P-T prot. - Lysis Buf.4 (4X) 130 mL	14RA
HTRF P-T prot. - Detect. Buf. 50 mL	33A
HTRF P-T prot. - Block. reag.(100X) 24 mL	26RA
HTRF EGFR total kit - 10K pts d2 antibody	12RA
HTRF EGFR total kit - 10K pts Eu Cryptate antibody	12RA
HTRF EGFR total kit - Ctrl lysate	09A

Results :

Incubation : Overnight at RT
Plate : 384 wells small volume white plate 20 µL
Reader : Envision
Fitting model : No curve

Assay characteristics :

	Found value	Specifications Rev 01
NEGATIF	397	/
Control Lysate	2292	/
S/B	5.8	>5
	/	/

HTRF EGFR total kit - 10K pts - Lot n° : 12G

☒ Accepted☐ Rejected

Quality control date : 2024-11-07

Supervisor's initials : PR

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. 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.