

HTRF® Research Reagents

HTRF EGFR p-pan kit - 500 pts

Product Number : 64HR1PEG Lot Number : 17E

Expiration date : 2025-07-01 Document version : 1

Reagents :	Lot n°
HTRF EGFR p-pan kit - Ctrl lysate	09A
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	24A
HTRF EGFR p-pan kit - 500 pts XL665 antibody	17B
HTRF EGFR p-pan kit - 500 pts Tb Cryptate antibody	08RA
HTRF EGFR- Lysis buffer (4X) - 2 mL	12A
HTRF EGFR Detect. Buf. - 2 mL	18A
HTRF EGFR-Stim. Buf. (5X) - 1.5 mL	15RA

Results :

Incubation : Overnight at RT  
 Plate : 384 wells small volume white plate 20 µL  
 Reader : Pherastar Plus  
 Fitting model : /

Assay characteristics :

	Found value	Specifications Rev 01
Negative	1124	/
Control Lysate	15398	/
S/B	13.7	>5
	/	/

HTRF EGFR p-pan kit - 500 pts - Lot n° : 17E

Accepted

Rejected

Quality control date : 2023-06-12	Supervisor's initials : CS
-----------------------------------	----------------------------

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.