

HTRF® Research Reagents

HTRF EGFR p-pan kit - 10K pts

Product Number : 64HR1PEH Lot Number : 16F

Expiration date : 2025-06-23 Document version : 1

Reagents :	Lot n°
HTRF EGFR p-pan kit - Ctrl lysate	08RA
HTRF P-T prot. - Block. reag.(100X) 2 mL	28A
HTRF EGFR p-pan kit - 10K pts XL665 antibody	16RA
HTRF EGFR p-pan kit - 10K pts Tb Cryptate antibody	07RA
HTRF EGFR- Lysis buffer (4X) - 130 mL	12RA
HTRF EGFR Detect. Buf. - 50 mL	19RA
HTRF EGFR-Stim. Buf. (5X) - 40 mL	15RA
HTRF EGFR p-pan kit - modop	8

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	1141	/
Control Lysate	9220	/
S/B	8.1	>5

HTRF EGFR p-pan kit - 10K pts - Lot n° : 16F

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.