

HTRF® Research Reagents

HTRF Adv.ERK p-T202/Y204 kit - 10K pts

Product Number : 64AERPEH Lot Number : 21B

Expiration date : 2026-04-03 Document version : 1

Reagents :	Lot n°
HTRF Adv.ERK p-T202/Y204 kit - 10K pts d2 antibody	21A
HTRF Adv.ERK p-T202/Y204 kit - 10K pts Eu Cryptate antibody	21A
HTRF Adv.ERK p-T202/Y204 kit - Ctrl lys.	20A
HTRF P-T prot. - Lysis Buf.1 (4X) 130 mL	34RA
HTRF P-T prot. - Block. reag.(100X) 2 mL	28B
HTRF P-T prot. - Block. reag.(100X) 2 mL	28B
HTRF P-T prot. - Detect. Buf. 50 mL	32RA

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	388	/
Control Lysate	2485	/
S/B	6,4	>5

HTRF Adv.ERK p-T202/Y204 kit - 10K pts - Lot n° : 21B

 Accepted

 Rejected

Quality control date : 2024-12-16	Supervisor's initials : ML
-----------------------------------	----------------------------

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.