

HTRF® Research Reagents

HTRF Adv.ERK p-T202/Y204 kit - 96 pts

Product Number : 64AERPET Lot Number : 17G

Expiration date : 2025-01-27 Document version : 1

Reagents :	Lot n°
PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND	2203
HTRF Adv.ERK p-T202/Y204 kit - Ctrl lys.	19A
HTRF Adv.ERK p-T202/Y204 kit - 96 pts d2 antibody	17RB
HTRF Adv.ERK p-T202/Y204 kit - 96 pts Eu Cryptate antibody	17RB
HTRF P-T prot. - Lysis Buf.1 (4X) 2 mL	35B
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	28RA
HTRF P-T prot. - Detect. Buf. 0.5 mL	23RB

Results :

Incubation :

Plate :

Reader : Pherastar Plus

Fitting model :

Assay characteristics :

	Found value	Specifications Rev 01
Negative	364	/
Control Lysate	2537	/
S/B	7	>5
	/	/

HTRF Adv.ERK p-T202/Y204 kit - 96 pts - Lot n° : 17G
 Accepted

 Rejected

Quality control date : 2024-03-14	Supervisor's initials : AC
-----------------------------------	----------------------------

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.