

## HTRF® Research Reagents

**HTRF ERK p-T202/Y204 kit - 500 pts**

Product Number : 64ERKPEG Lot Number : 19M

Expiration date : 2026-01-11 Document version :

Reagents :	Lot n°
HTRF P-T prot. - Lysis Buf.1 (4X) 8 mL	35AR
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	29A
HTRF P-T prot. - Detect. Buf. 4 mL	32A
HTRF ERK p-T202/Y204 kit - Ctrl lysate	15RA
HTRF ERK p-T202/Y204 kit - 500 pts d2 antibody	19B
HTRF ERK p-T202/Y204 kit - 500 pts Eu Cryptate antibody	19B

**Results :**

 Incubation : 2h at RT  
 Plate : 384 wells small volume white plate 20 µL  
 Reader : EnVision  
 Fitting model : /

**Assay characteristics :**

	Found value	Specifications Rev 01
Negative	527	/
Control Lysate	2991	/
S/B	5.7	>5

HTRF ERK p-T202/Y204 kit - 500 pts - Lot n° : 19M

 Accepted

 Rejected

Quality control date : 2024-11-26	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.