

HTRF® Research Reagents

HTRF P38 p-T180/Y182 kit - 96 pts
Product Number : 64P38PET **Lot Number :** 01ZA

Expiration date : 09/03/2025 **Document version :** 1

Reagents :	Lot n°
PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND	2203
HTRF P38 p-T180/Y182 kit - Ctrl lysate	12RA
HTRF P38 p-T180/Y182 kit - 96 pts d2 antibody	01RA
HTRF P38 p-T180/Y182 kit - 96 pts Eu Cryptate antibody	01RA
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 : 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	538	/
Control Lysate	6796	/
S/B	12,6	>5
	/	/

HTRF P38 p-T180/Y182 kit - 96 pts - Lot n° : 01ZA
 Accepted

 Rejected

Quality control date :	2023-01-17	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.