

HTRF® Research Reagents

HTRF AMPK p-T172 kit - 96 pts

Product Number : 64MPKPET Lot Number : 01X

Expiration date : 2024-12-17 Document version : 1

Reagents :	Lot n°
PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND	2203
PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND	2203
HTRF P-T prot. - Lysis Buf.1 (4X) 2 mL	33A
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	28A
HTRF P-T prot. - Detect. Buf. 0.5 mL	23RB
HTRF AMPK p-T172 kit - Ctrl lysate	12RA
HTRF AMPK p-T172 kit - 96 pts modop	5
HTRF AMPK p-T172 kit - 96 pts d2 antibody	11RC
HTRF AMPK p-T172 kit - 96 pts Eu Cryptate antibody	11RC

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	428	/
Control Lysate	3580	/
S/B	8.4	>5
	/	/

HTRF AMPK p-T172 kit - 96 pts - Lot n° : 01X

 Accepted

 Rejected

Quality control date : 2022-10-12	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.