

HTRF® Research Reagents

HTRF AMPK total kit - 500 pts

Product Number : 63ADK060PEG Lot Number : 11B

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

Reagents :	Lot n°
HTRF AMPK total kit - 500 pts d2 antibody	11B
HTRF AMPK total kit - 500 pts Eu Cryptate antibody	11B
HTRF AMPK total kit - Ctrl lysate	09A
HTRF P-T prot. - Lysis Buf.1 (4X) 2 mL	35AR
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	28RA
HTRF P-T prot. - Detect. Buf. 2 mL	32B

Results :

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

Assay characteristics :

	Found value	Specifications Rev 01
Negative	417	/
Control Lysate	3695	/
S/B	8.9	>5
	/	/

HTRF AMPK total kit - 500 pts - Lot n° : 11B

☒ Accepted

☐ Rejected

Quality control date : 2024-02-16 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.