

HTRF® Research Reagents

## HTRF AKT2 p-S473 kit - 10K pts

Product Number : 63ADK080PEH Lot Number : 11C

Expiration date : 06/03/2025 Document version : 1

Reagents :	Lot n°
HTRF AKT2 p-S473 kit - 10K pts d2 antibody	11A
HTRF AKT2 p-S473 kit - 10K pts Eu Cryptate antibody	11A
HTRF AKT2 p-S473 kit - Ctrl lysate	12AR
HTRF AKT2 p-S473 kit - modop	8
HTRF P-T prot. - Lysis Buf.1 (4X) 130 mL	34RA
HTRF P-T prot. - Block. reag.(100X) 2 mL	28A
HTRF P-T prot. - Detect. Buf. 50 mL	32RA

**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	432	/
Control Lysate	3137	/
S/B	7,3	>5
	/	/

HTRF AKT2 p-S473 kit - 10K pts - Lot n° : 11C

 Accepted

 Rejected

Quality control date : 2023-09-06	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.