

## HTRF® Research Reagents

**HTRF AKT p-S473 kit - 96 pts**

Product Number : 64AKSPET Lot Number : 19G

Expiration date : 2025-01-13 Document version : 1

Reagents :	Lot n°
PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND	2203
HTRF AKT p-S473 kit - Ctrl lysate	19RA
HTRF AKT p-S473 kit - 96 pts modop	5
HTRF AKT p-S473 kit - 96 pts d2 antibody	19RA
HTRF AKT p-S473 kit - 96 pts Eu Cryptate antibody	19RA
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

**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	299	/
Control Lysate	5889	/
S/B	19.7	>5
	/	/

HTRF AKT p-S473 kit - 96 pts - Lot n° : 19G

 Accepted

 Rejected

Quality control date : 2023-01-09	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.