

HTRF® Research Reagents

HTRF AKT p-S473 kit - 500 pts

Product Number : 64AKSPEG Lot Number : 19K

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

Reagents :	Lot n°
HTRF AKT p-S473 kit - 500 pts d2 antibody	19RA
HTRF AKT p-S473 kit - 500 pts Eu Cryptate antibody	19RA
HTRF AKT p-S473 kit - Ctrl lysate	19RA
HTRF AKT p-S473 kit - modop	7
HTRF P-T prot. - Lysis Buf.1 (4X) 2 mL	34A
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	28A
HTRF P-T prot. - Detect. Buf. 2 mL	32A

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	329	/
Control Lysate	5367	/
S/B	16.3	>5
	/	/

HTRF AKT p-S473 kit - 500 pts - Lot n° : 19K

Accepted

Rejected

Quality control date : 2023-01-09	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.