

HTRF® Research Reagents

## HTRF AKT p-S473 kit - 500 pts

Product Number : 64AKSPEG Lot Number : 20B

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

Reagents :	Lot n°
HTRF AKT p-S473 kit - 500 pts d2 antibody	20A
HTRF AKT p-S473 kit - 500 pts Eu Cryptate antibody	20A
HTRF AKT p-S473 kit - Ctrl lysate	20AR
HTRF P-T prot. - Lysis Buf.1 (4X) 2 mL	35B
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	28B
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	307	/
Control Lysate	5133	/
S/B	16.7	>5
	/	/

HTRF AKT p-S473 kit - 500 pts - Lot n° : 20B

 Accepted

 Rejected

Quality control date : 2024-04-24 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.