

HTRF® Research Reagents

HTRF AKT p-S473 kit - 10K pts

Product Number : 64AKSPEH Lot Number : 21C

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

Reagents :	Lot n°
HTRF AKT p-S473 kit - Ctrl lysate	20AR
HTRF AKT p-S473 kit - 10K pts d2 antibody	21A
HTRF AKT p-S473 kit - 10K pts Eu Cryptate antibody	21A
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 : Envision
 Fitting model : /

Assay characteristics :

	Found value	Specifications Rev 01
Negative	300	/
Control Lysate	3042	/
S/B	10.1	>5
	/	/

HTRF AKT p-S473 kit - 10K pts - Lot n° : 21C

 Accepted

 Rejected

Quality control date : 2023-10-17	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.