

HTRF® Research Reagents

HTRF YAP p-S127 kit - 500 pts

Product Number : 64YAPPEG Lot Number : 11A

Expiration date : 2026-01-07 Document version : 1

Reagents :	Lot n°
HTRF P-T prot. - Lysis Buf.4 (4X) 2 mL	14A
HTRF P-T prot. - Block. reagent.(100X) 0.3 mL	28A
HTRF P-T prot. - Detect. Buf. 2 mL	32A
HTRF YAP p-S127 kit - 500 pts d2 antibody	11B
HTRF YAP p-S127 kit - 500 pts Eu Cryptate antibody	11B
HTRF YAP p-S127 kit - Ctrl lysate	10RA
HTRF YAP p-S127 kit - modop	5

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	374	/
Control Lysate	3508	/
S/B	9.4	>5
	/	/

HTRF YAP p-S127 kit - 500 pts - Lot n° : 11A

☒ Accepted☐ Rejected

Quality control date : 2023-03-22	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.