

HTRF® Research Reagents

HTRF (h) YES Total Kit - 500 pts

Product Number : 64YESTPEG Lot Number : 01A

Expiration date : 2027-04-14 Document version : 1

Reagents :	Lot n°
HTRF (h) YES Total Kit - 500 pts d2 antibody	01A
HTRF (h) YES Total Kit - 500 pts Eu Cryptate antibody	01A
HTRF (h) YES Total Kit - Ctrl Lysate	01A
HTRF P-T prot. - Lysis Buf.4 (4X) 2 mL	14B
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	29A
HTRF P-T prot. - Detect. Buf. 2 mL	33A

Results :

Incubation : 4h at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Envision
 Fitting model : No curve

Assay characteristics :

	Found value	Specifications Rev 01
NEGATIF	411	/
Control Lysate	3943	/
S/B	9.6	>5
	/	/

HTRF (h) YES Total Kit - 500 pts - Lot n° : 01A

☒ Accepted

☐ Rejected

Quality control date : 2025-01-13 Supervisor's initials : PR

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.