

HTRF® Research Reagents

HTRF (h) SMARCA2 Total Kit - 500 pts

Product Number : 64SMARC2TPEG Lot Number : 02b

Expiration date : 2025-08-04 Document version : 1

Reagents :		Lot n°
HTRF (h) SMARCA2 Total Kit - 500 pts d2 antibody		02A
HTRF (h) SMARCA2 Total Kit - 500 pts Eu Cryptate antibody		02A
HTRF (h) SMARCA2 Total Kit - Ctrl Lys		01RA
HTRF P-T prot. - Lysis Buf.1 (4X) 2 mL		35B
HTRF P-T prot. - Block. reag.(100X) 0.3 mL		28RA
HTRF P-T prot. - Detect. Buf. 2 mL		32B

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	503	
Control Lysate	5293	/
S/B	10.5	>5
	/	/

HTRF (h) SMARCA2 Total Kit - 500 pts - Lot n° : 02B

☒ Accepted☐ Rejected

Quality control date : 2023-02-16 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.