

HTRF® Research Reagents
HTRF RB p-S807/811 kit - 500 pts

Product Number : 64RBS807PEG Lot Number : 14B

Expiration date : 2027-06-06 Document version : 1

Reagents :	Lot n°
HTRF RB p-S807/811 kit - 500 pts d2 antibody	14A
HTRF RB p-S807/811 kit - 500 pts Eu Cryptate antibody	14A
HTRF RB p-S807/811 kit - Ctrl lysate	09A
HTRF P-T prot. - Lysis Buf.2 (4X) 2 mL	13AR
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	29A
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 : /

Assay characteristics :

	Found value	Specifications Rev 01
Negative	339	/
Control Lysate	4088	/
S/B	11	>5

HTRF RB p-S807/811 kit - 500 pts - Lot n° : 14B

☒ Accepted

☐ Rejected

Quality control date : 2025-01-31	Supervisor's initials : ML
-----------------------------------	----------------------------

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.