

HTRF® Research Reagents

## HTRF RB total kit - 10K pts

Product Number : 64RBTPPEH Lot Number : 07A

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

Reagents :	Lot n°
HTRF RB total kit - 10K pts d2 antibody	07A
HTRF RB total kit - 10K pts Eu Cryptate antibody	07A
HTRF RB total kit - Ctrl lysate	05A
HTRF P-T prot. - Lysis Buf.2 (4X) 130 mL	13AR
HTRF P-T prot. - Block. reag.(100X) 2 mL	28B
HTRF P-T prot. - Detect. Buf. 50 mL	32RA

**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	363	/
Control Lysate	2923	/
S/B	8	>5
	/	/

HTRF RB total kit - 10K pts - Lot n° : 07A

 Accepted

 Rejected

Quality control date : 2024-05-30 Supervisor's initials : AC

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.