

HTRF® Research Reagents

HTRF RB total kit - 500 pts

 Product Number : **64RBTPEG** Lot Number : **06C**

 Expiration date : **2025-02-24** Document version : **1**

Reagents :	Lot n°
HTRF RB total kit - 500 pts d2 antibody	06AR
HTRF RB total kit - 500 pts Eu Cryptate antibody	06AR
HTRF RB total kit - Ctrl lysate	04RA
HTRF P-T prot. - Lysis Buf.2 (4X) 2 mL	12RA
HTRF P-T prot. - Block. reag.(100X) 0.3 mL	26RA
HTRF P-T prot. - Detect. Buf. 2 mL	32A

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	383	/
Control Lysate	3952	/
S/B	10,3	>5
	/	/

HTRF RB total kit - 500 pts - Lot n° : 06C
 Accepted

 Rejected

Quality control date : 2023-11-07	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.