

HTRF® Research Reagents

HTRF B-arrestin 1 total kit - 1K pts

Product Number : 64BAR1TPEB Lot Number : 01C

Expiration date : 2025-01-15 Document version : 1

Reagents :	Lot n°
HTRF Lysis Buffer 4 - 16 mL	12RA
HTRF B-arrestin 1 total kit - 1K pts d2 antibody	01RA
HTRF B-arrestin 1 total kit - 1K pts Eu Cryptate antibody	01RA
HTRF B-arrestin 1 total kit - Ctrl lys.	01RA
HTRF P-T prot. - Detec. 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
Control Lysate	4476	> 4300
Ratio negative	389	380-449
S/B	11,5	/
	/	/

HTRF B-arrestin 1 total kit - 1K pts - Lot n° : 01C

Accepted

Rejected

Quality control date : 2024-04-05	Supervisor's initials : MP
-----------------------------------	----------------------------

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.