

HTRF® Research Reagents
HTRF B-arrestin 1 total kit - 1K pts

Product Number : 64BAR1TPEB Lot Number : 01D

Expiration date : 2025-11-25 Document version : 1

Reagents :	Lot n°
HTRF Lysis Buffer 4 - 16 mL	14RA
HTRF B-arrestin 1 total kit - 1K pts d2 antibody	01RA
HTRF B-arrestin 1 total kit - 1K ptsEu Cryptate antibody	01RA
HTRF B-arrestin 1 total kit - Ctrl lys.	01RA
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 : No curve

Assay characteristics :

	Found value	Specifications Rev 01
Negative	4674	> 4300
Control Lysate	418	380-449
S/B	11	

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

☒ Accepted

☐ Rejected

Quality control date : 2026-12-24	Supervisor's initials : FMV
-----------------------------------	-----------------------------

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.