

HTRF® Research Reagents

HTRF B-arr2 recruitment kit - 1K pts

Product Number : 62BDBAR2PEB Lot Number : 06E

Expiration date : 2025-09-01 Document version : 1

Reagents :	Lot n°
HTRF Stimulation Buf. 4 (5X) - 20 mL	15RB
HTRF Stabilization Buf. 1 - 30 mL	02RA
HTRF Wash Buffer 1 (10X) - 30 mL	02RA
HTRF Detect. Buf. 14 (10X) - 10 mL	02RB
HTRF B-arr2 recruitment kit - 1K pts d2 antibody	06RA
HTRF B-arr2 recruitment kit - 1K pts Eu Cryptate antibody	06RA
HTRF B-arr2 recruitment kit - modop	2

Results :

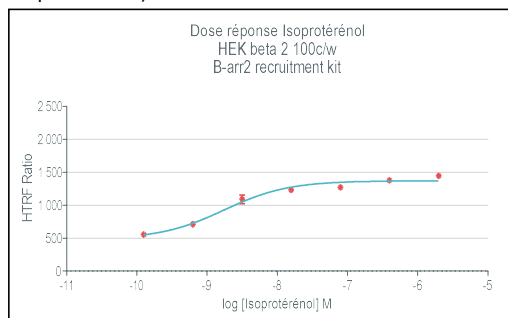
Incubation : overnight dans étuve thermostatée (20°C)

Plate : Plaque 96 cultureplate

Reader : Pherastar Plus

Fitting model : log(agonist)vs.response. Variable slope (four parameters)

Agonist	Conc. (Mf)	ratio
no ago	basal	526
agonist 1	1,3E-01	555
agonist 2	6,4E-01	711
agonist 3	3,2E+00	1091
agonist 4	1,6E+01	1233
agonist 5	8,0E+01	1272
agonist 6	4,0E+02	1379
agonist 7	2,0E+03	1447



Assay characteristics :

	Found value	Specifications Rev 01
EC50	17 **	20 - 80
S/B	2,7	1.4 - 3.6
	/	/
	/	/

** Batch validated with EC50 a little bit lower than spécification

HTRF B-arr2 recruitment kit - 1K pts - Lot n° : 06E

Accepted

Rejected

Quality control date : 2023-09-14 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.