

HTRF® Research Reagents

HTRF B-arr2 recruitment kit - 1K pts

Product Number : 62BDBAR2PEB Lot Number : 07A

Expiration date : 2026-03-17 Document version : 1

Reagents :	Lot n°
HTRF Stimulation Buf. 4 (5X) - 20 mL	18RA
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	07A
HTRF B-arr2 recruitment kit - 1K pts Eu Cryptate antibody	07A

Results :

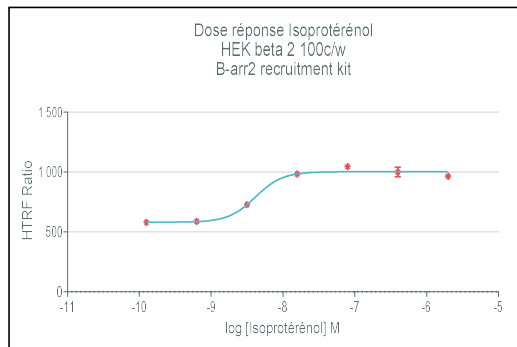
Incubation : overnight dans étuve thermostatée (20°C)

Plate : Plaque 96 cultureplate

Reader : Envision

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

Agonist	Conc. (Mf)	ratio
no ago	basal	586
agonist 1	1,3E-01	580
agonist 2	6,4E-01	588
agonist 3	3,2E+00	728
agonist 4	1,6E+01	984
agonist 5	8,0E+01	1045
agonist 6	4,0E+02	1001
agonist 7	2,0E+03	965



Assay characteristics :

	Found value	Specifications Rev 01
EC50	4,1	20 - 80
S/B	1,6	1.4 - 3.6
	/	/
	/	/

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

HTRF B-arr2 recruitment kit - 1K pts - Lot n° : 07A

Accepted

Rejected

Quality control date : 2024-02-28	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.