

HTRF® Research Reagents

HTRF B-arr2 recruitment kit - 1K pts

Product Number : 62BDBAR2PEB Lot Number : 08B

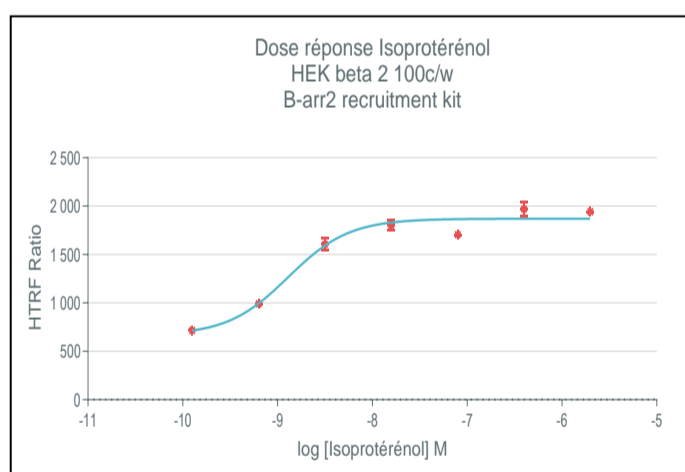
Expiration date : 2026-03-14 Document version : 1

Reagents :	Lot n°
HTRF Stimulation Buf. 4 (5X) - 20 mL	18RA
HTRF Stabilization Buf. 1 - 30 mL	03A
HTRF Wash Buffer 1 (10X) - 30 mL	02RA
HTRF Detect. Buf. 14 (10X) - 10 mL	02C
HTRF B-arr2 recruitment kit - 1K pts d2 antibody	08A
HTRF B-arr2 recruitment kit - 1K pts Eu Cryptate antibody	08A

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	581
agonist 1	1,3E-01	715
agonist 2	6,4E-01	991
agonist 3	3,2E+00	1606
agonist 4	1,6E+01	1806
agonist 5	8,0E+01	1706
agonist 6	4,0E+02	1969
agonist 7	2,0E+03	1943



Assay characteristics :

	Found value	Specifications Rev 01
EC50	13	20 - 80
S/B	3,3	1.4 - 3.6
	/	/
	/	/

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

HTRF B-arr2 recruitment kit - 1K pts - Lot n° : 08B

Accepted

Rejected

Quality control date : 2024-11-27	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.