

HTRF® Research Reagents

HTRF IL23/IL12Rb1 Bind. kit - 500 pts

Product Number : 64BDPIL12PEG Lot Number : 02B

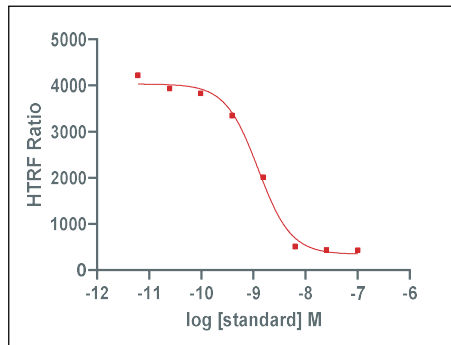
Expiration date : 2025-11-19 Document version : 1

Reagents :		Lot n°
HTRF PPI-Europium Detect. Buf. - 20 mL		07RA
HTRF IL23/IL12Rb1 Bind. kit - 500 pts Tag2-IL12Rb1		02RA
HTRF IL23/IL12Rb1 Bind. kit - Std		02RA
HTRF IL23/IL12Rb1 Bind. kit - 500 pts Eu Cryptate antibody		02RA
HTRF IL23/IL12Rb1 Bind. kit - 500 pts Tag1-IL23		01RA
HTRF IL23/IL12Rb1 Bind. kit - 500 pts XL665 reagent		02RA

Results :

Incubation : 2h at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model : log(inhibitor) vs. response -- Variable slope (four parameters)

std	Conc. (pM)	RATIO
Std 0	0	4223
Std 1	24,4	3937
Std 2	97,7	3830
Std 3	390	3349
Std 4	1563	2014
Std 5	6250	509
Std 6	25000	437
Std 7	100000	430



Assay characteristics :

	Found value	Specifications Rev 01
negatif	450	
S/B	9	8 -15
IC50	1,25	0.7 -2 nM

HTRF 1L23/IL12Rb1 Bind, kit - 500 pts Lot 02B

Accepted

Rejected

Quality control date : 2024-04-09 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 TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.