

HTRF® Research Reagents

HTRF IL23/IL23R Bind. kit - 500 pts

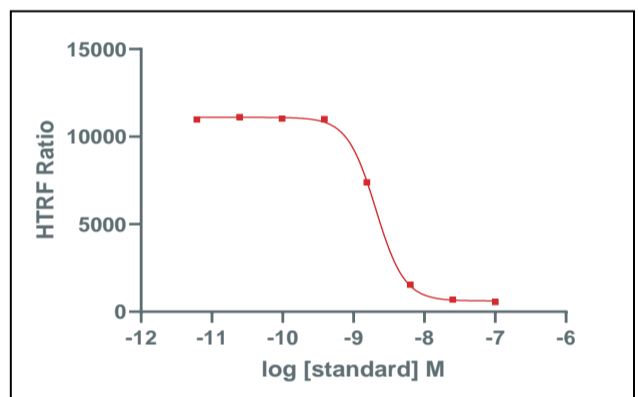
Product Number : 64BDPIL23PEG Lot Number : 03K

Expiration date : 2025-07-28 Document version : 1

Reagents :	Lot n°
HTRF PPI-Europium Detect. Buf. - 20 mL	07RA
HTRF IL23/IL23R Bind. kit - 500 pts Tag2-IL23R	05RA
HTRF IL23/IL23R Bind. kit - Std	02RA
HTRF IL23/IL23R Bind. kit - 500 pts d2 reagent	04A
HTRF IL23/IL23R Bind. kit - 500 pts Eu Cryptate antibody	04A
HTRF IL23/IL23R Bind. kit - 500 pts Tag1-IL23	05RA
HTRF IL23/IL23R Bind. kit - modop	1

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

std	Conc. (pM)	RATIO
Std 0	0	10 982
Std 1	24,4	11 124
Std 2	97,7	11 034
Std 3	390	11 009
Std 4	1563	7 390
Std 5	6250	1 543
Std 6	25000	693
Std 7	100000	554



Assay characteristics :

	Found value	Specifications Rev 01
negatif	392	/
S/B	28	15-30
IC50	2,06	1.5 - 3.4 nM

HTRF IL23/IL23R Bind. kit - 500 pts - Lot n° : 03K

Accepted

Rejected

Quality control date : 2023-05-25	Supervisor's initials : PR
-----------------------------------	----------------------------

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.