

HTRF® Research Reagents

HTRF IL23/IL23R Bind. kit - 500 pts

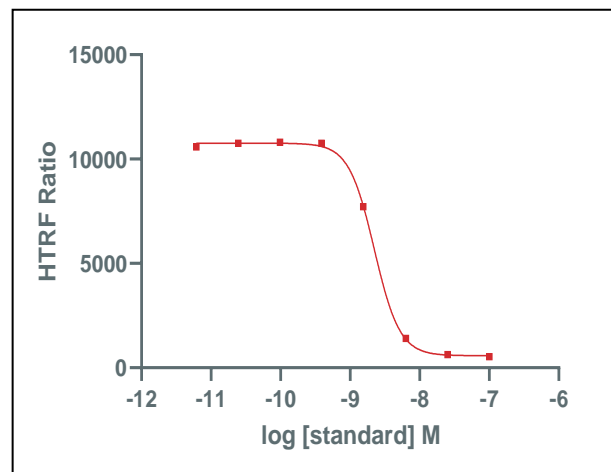
Product Number : 64BDPIL23PEG Lot Number : 03H

Expiration date : 2025-07-25 Document version : 1

Reagents :	Lot n°
HTRF PPI-Europium Detect. Buf. - 20 mL	07RA
HTRF IL23/IL23R Bind. kit - 500 pts Tag2-IL23R	03RA
HTRF IL23/IL23R Bind. kit - Std	02RA
HTRF IL23/IL23R Bind. kit - 500 pts d2 reagent	03RA
HTRF IL23/IL23R Bind. kit - 500 pts Eu Cryptate antibody	03RA
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	10578
Std 1	24,4	10754
Std 2	97,7	10809
Std 3	390	10754
Std 4	1563	7719
Std 5	6250	1404
Std 6	25000	632
Std 7	100000	533



Assay characteristics :

	Found value	Specifications Rev 01
negatif	390	/
S/B	27	15-30
IC50	2,2	1.5 - 3.4 nM

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

Accepted

Rejected

Quality control date : 2023-09-13	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.