

HTRF® Research Reagents

HTRF Human IL23 kit - 500 pts

Product Number : 62HIL23PEG Lot Number : 06B

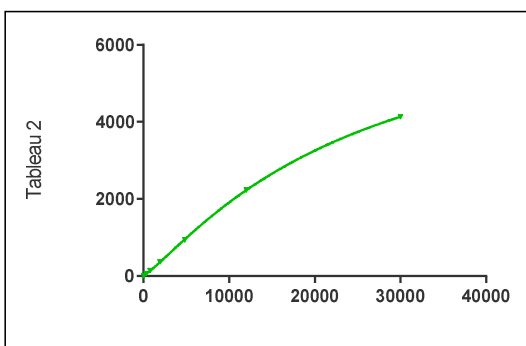
Expiration date : 2025-08-17 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 1.5 mL	30AR
HTRF Human IL23 kit - Std	04RA
HTRF Human IL23 kit - 500 pts d2 antibody	06AR
HTRF Human IL23 kit - 500 pts Eu Cryptate antibody	06AR
HTRF Human IL23 kit - modop	7

Results :

Incubation : **3h at 4°C**
 Plate : **384 wells small volume white plate 20 µL**
 Reader : **Pherastar Plus**
 Fitting model : **4 Parameter Logistic (4PL 1/y2)* model**

IL23	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	123	24%	16-38
Std 2	307	61%	46-93
Std 3	768	142%	122-183
Std 4	1920	367%	317-457
Std 5	4800	936%	759-1146
Std 6	12000	2233%	1828-2383
Std 7	30000	4133%	3528-4597



Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	20 pg/ml	≤ 31 pg/ml

HTRF Human IL23 kit - 500 pts - Lot n° : 06B

Accepted

Rejected

Quality control date :	2023-11-21	Supervisor's initials :	PL
------------------------	------------	-------------------------	----

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.