

HTRF® Research Reagents

HTRF (h) IL12/23 p40 Kit - 500 pts

Product Number : 62HIL12P40PEG Lot Number : 03A

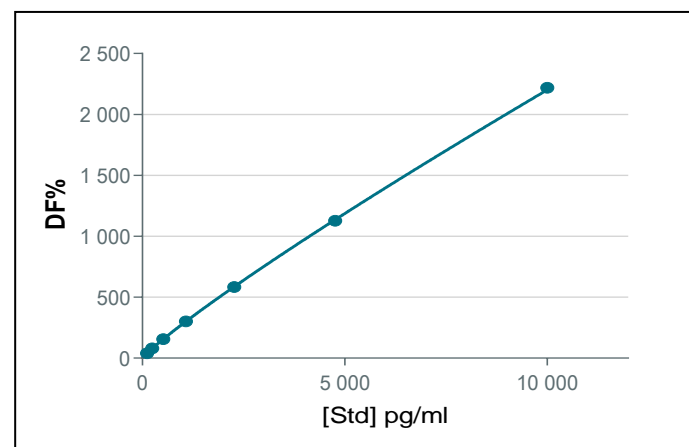
Expiration date : 2025-10-28 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21B
HTRF Detect. Buf. 3 - 1.5 mL	30B
HTRF (h) IL12/23 p40 Kit - Std	01RA
HTRF (h) IL12/23 p40 Kit - 500 pts d2 antibody	03A
HTRF (h) IL12/23 p40 Kit - 500 pts Eu Cryptate antibody	03A
HTRF (h) IL12/23 p40 Kit - modop	1

Results :

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

h IL12P40 Kit	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	117	40%	28 --- 40
Std 2	245	80%	52 --- 80
Std 3	514	155%	118 --- 159
Std 4	1080	301%	235 --- 310
Std 5	2268	584%	541 --- 607
Std 6	4762	1127%	1078 --- 1262
Std 7	10000	2219%	2162 --- 2528



Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	11,6 pg/ml	< 117 pg/ml

HTRF (h) IL12/23 p40 Kit - 500 pts - Lot n° : 03A

Accepted

Rejected

Quality control date : 2024-08-23	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.