

HTRF® Research Reagents

HTRF HP Human IL6 Kit - 500 pts

Product Number : 62HIL6V2PEG Lot Number : 02A

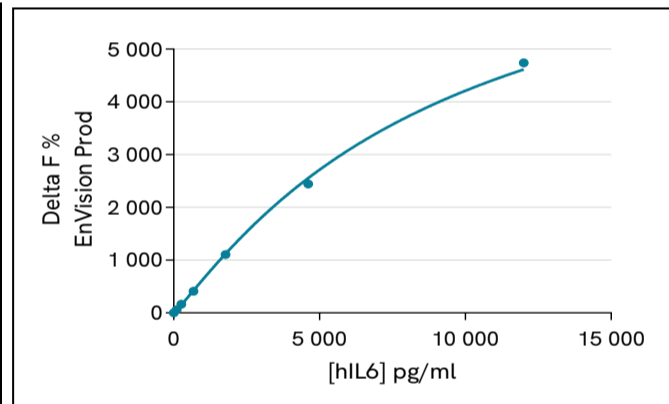
Expiration date : 2027-05-23 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21B
HTRF Detect. Buf. 3 - 1.5 mL	30B
HTRF HP Human IL6 Kit - Std	01A
HTRF HP Human IL6 Kit - 500 pts d2 antibody	02A
HTRF HP Human IL6 Kit - 500 pts Eu Cryptate antibody	02A

Results :

Incubation : 2h at RT  
 Plate : 384 wells small volume white plate 20 µL  
 Reader : Envision Lamp  
 Fitting model : 4 Parameter Logistic (4PL 1/y2)\* model

hIL6	Conc. (pg/ml)	DF %	DF % Accepted Range
Std 1	38,8	17%	12 - 43
Std 2	101	59%	35 - 90
Std 3	262,6	164%	111 - 211
Std 4	682,7	410%	312 - 473
Std 5	1775,1	1106%	791 - 1186
Std 6	4615,5	2444%	1897 - 2845
Std 7	12000	4740%	4031 - 6047



Assay characteristics :

	Found value	Specifications Rev 01
Limite de detection	8,9 pg/ml	< 37 pg/ml

HTRF HP Human IL6 Kit - 500 pts - Lot n° : 02A

Accepted

Rejected

Quality control date : 2025-01-23	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.