

HTRF® Research Reagents

HTRF HP Human IL6 Kit - 10K pts

Product Number : 62HIL6V2PEH Lot Number : 01A

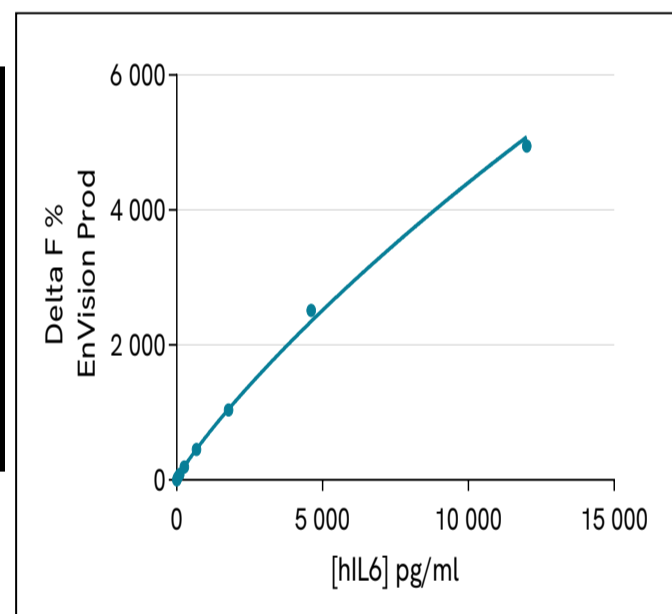
Expiration date : 2027-04-05 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 10 mL	21RA
HTRF Detect. Buf. 3 - 50 mL	29RA
HTRF HP Human IL6 Kit - 10K pts d2 antibody	01A
HTRF HP Human IL6 Kit - 10K pts Eu Cryptate antibody	01A
HTRF HP Human IL6 Kit - Std	01A

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	39	12 - 43
Std 2	101	83	35 - 90
Std 3	262,6	192	111 - 211
Std 4	682,7	454	312 - 473
Std 5	1775,1	1035	791 - 1186
Std 6	4615,5	2510	1897 - 2845
Std 7	12000	4948	4031 - 6047



Assay characteristics :

	Found value	Specifications Rev 01
LoD	19,2 pg/ml	< 37 pg/ml

HTRF HP Human IL6 Kit - 10K pts - Lot n° : 01A

Accepted

Rejected

Quality control date : 2024-07-18	Supervisor's initials : FL
-----------------------------------	----------------------------

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.