

HTRF® Research Reagents

HTRF Human IL6 kit - 500 pts

Product Number : 62HIL06PEG Lot Number : 13A

Expiration date : 2027-07-26 Document version : 1

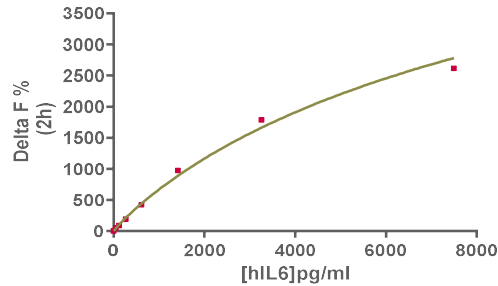
Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21B
HTRF Detect. Buf. 3 - 1.5 mL	30B
HTRF Human IL6 kit - Std	07A
HTRF Human IL6 kit - 500 pts Eu Cryptate antibody	13A
HTRF Human IL6 kit - 500 pts XL665 antibody	13A

Results :

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

Assay characteristics :

hIL6	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	50.7	53%	25-57
Std 2	117	90%	60-118
Std 3	268	195%	128-251
Std 4	616	421%	300-495
Std 5	1417	976%	632-1034
Std 6	3260	1787%	1255-2095
Std 7	7500	2614%	1875-3131



Assay characteristics :

	Found value	Specifications Rev 02
Detection limit	3.9 pg/ml	< 50.7 pg/ml

HTRF Human IL6 kit - 500 pts - Lot n° : 13A

Accepted

Rejected

Quality control date : 2024-08-05	Supervisor's initials : AC
-----------------------------------	----------------------------

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.