

HTRF® Research Reagents

HTRF Human IL6 kit - 500 pts

Product Number : 62HIL06PEG Lot Number : 09E

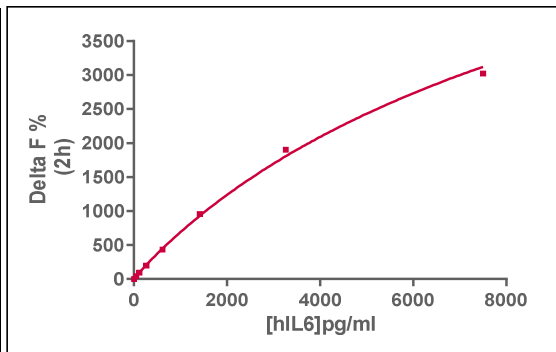
Expiration date : 2025-08-22 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 1.5 mL	30AR
HTRF Human IL6 kit - Std	06RA
HTRF Human IL6 kit - 500 pts Eu Cryptate antibody	09RA
HTRF Human IL6 kit - 500 pts XL665 antibody	09RA
HTRF Human IL6 kit - modop	7

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

hIL6	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	50.7	39%	25-57
Std 2	117	99%	60-118
Std 3	268	209%	128-251
Std 4	616	451%	300-495
Std 5	1417	973%	632-1034
Std 6	3260	1943%	1255-2095
Std 7	7500	3099%	1875-3131



Assay characteristics :

	Found value	Specifications Rev 02
Detection limit	5 pg/ml	< 50.7 pg/ml

HTRF Human IL6 kit - 500 pts - Lot n° : 09E

Accepted

Rejected

Quality control date :	2023-11-17	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.