

HTRF® Research Reagents

HTRF Human IL6 kit - 96 pts

Product Number : 62HIL06PET Lot Number : 12A

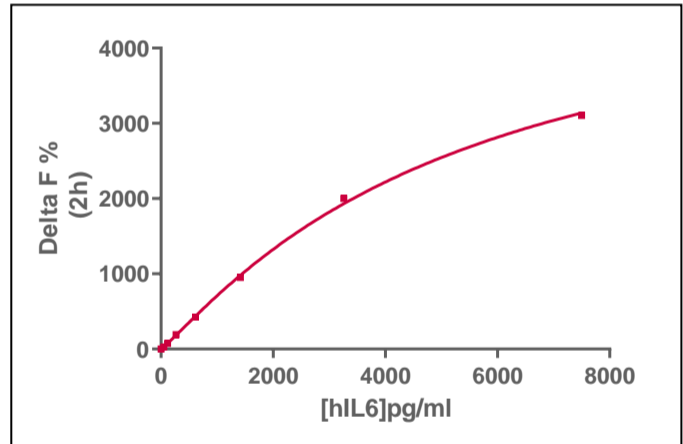
Expiration date : 2025-08-22 Document version : 1

Reagents :	Lot n°
PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND	2413
PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND	2413
HTRF Diluent 5 (5X) - 2 mL	21B
HTRF Detect. Buf. 3 - 0.5 mL	30AR
HTRF Human IL6 kit - Std	06RA
HTRF Human IL6 kit - 96 pts Eu Cryptate antibody	11A
HTRF Human IL6 kit - 96 pts XL665 antibody	12A

Results :

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

hIL6	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	50,7	29%	25-57
Std 2	117	75%	60-118
Std 3	268	191%	128-251
Std 4	616	424%	300-495
Std 5	1417	951%	632-1034
Std 6	3260	2001%	1255-2095
Std 7	7500	3105%	1875-3131



Assay characteristics :

	Found value	Specifications Rev 03
Detection limit	pg/ml	≤ 50.7 pg/ml

HTRF Human IL6 kit - 96 pts - Lot n° : 12A

Accepted

Rejected

Quality control date : 2024-07-02	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.