

HTRF® Research Reagents

HTRF HP Human TNF alpha Kit - 10K pts

Product Number : 62HTNFAV2PEH Lot Number : 01A

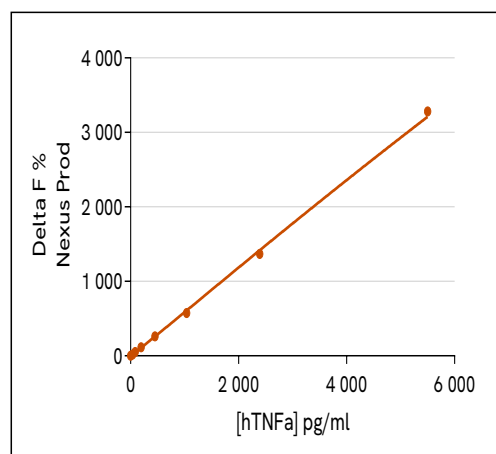
Expiration date : 2027-06-16 Document version : 1

Reagents :	Lot n°
HTRF Diluent 1 - 20 mL	20RA
HTRF Detect. Buf. 3 - 1.5 mL	29RA
HTRF HP Human TNF alpha Kit - Std	01A
HTRF HP Human TNF alpha Kit - 10K pts d2 antibody	01A
HTRF HP Human TNF alpha Kit - 10K pts Eu Cryptate antibody	01A

Results :

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

[hTNFa] pg/ml	Conc	DF%	DF % Accepted Range
Std 0	negatif (Std0)	0	0
Std 1	37,2	17	13 - 27
Std 2	85,5	53	42 - 62
Std 3	196,5	116	87 - 130
Std 4	452	261	191 - 287
Std 5	1039,7	575	404 - 606
Std 6	2391,3	1368	955 - 1433
Std 7	5500	3282	2362 - 3543



Assay characteristics :

	Found value	Specifications Rev 01
Limite de détection	8,5 pg/ml	< 37,2 pg/ml

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

☒ Accepted☐ Rejected

Quality control date : 2025-06-27 Supervisor's initials : MP

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.