

HTRF® Research Reagents

HTRF Human TNF alpha kit - 500 pts

Product Number : 62HTNFAPEG Lot Number : 12A

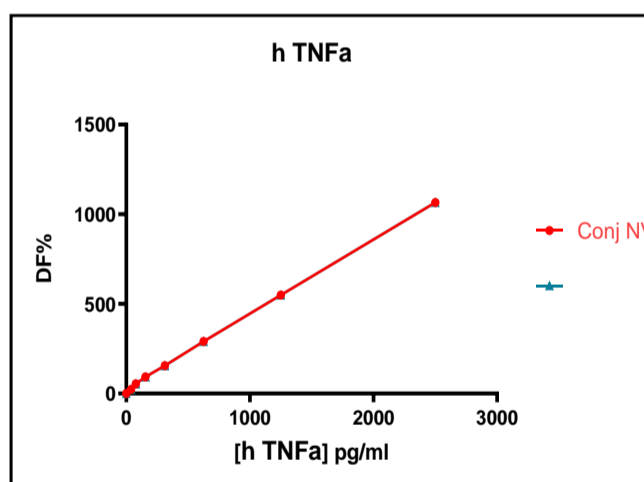
Expiration date : 2025-09-07 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21B
HTRF Detect. Buf. 3 - 1.5 mL	30B
HTRF Human TNF alpha kit - Std	05RA
HTRF Human TNF alpha kit - 500 pts d2 antibody	12A
HTRF Human TNF alpha kit - 500 pts Eu Cryptate 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

hTNF alpha	Conc. (pg/ml)	DF %	DF Accepted Range Revision 2
Std 1	39	23%	13-33
Std 2	78	54%	34-61
Std 3	156	93%	70-119
Std 4	312	155%	120-216
Std 5	625	290%	220-359
Std 6	1250	549%	400-645
Std 7	2500	1065%	730-1100



Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	6.8pg/ml	<39pg/ml

HTRF Human TNF alpha kit - 500 pts - Lot n° : 12A

Accepted

Rejected

Quality control date : 2024-10-07	Supervisor's initials : ML
-----------------------------------	----------------------------

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