

HTRF® Research Reagents

HTRF Human TNF alpha kit - 500 pts

Product Number : 62HTNFAPEG Lot Number : 09J

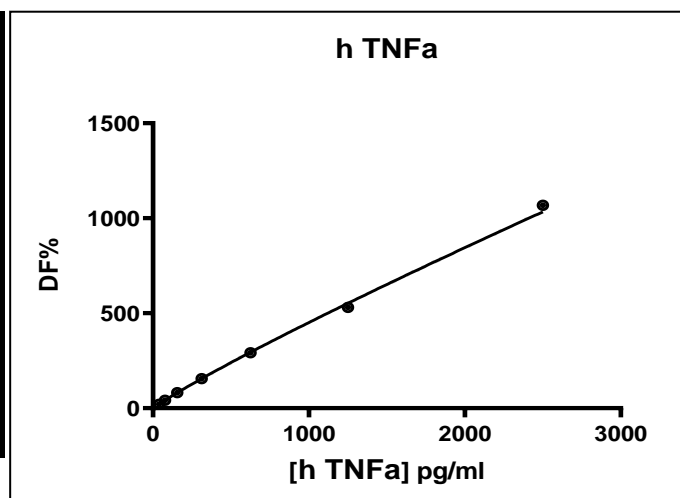
Expiration date : 2025-12-15 Document version : 1

Reagents :		Lot n°
HTRF Diluent 5 (5X) - 2 mL		21B
HTRF Detect. Buf. 3 - 1.5 mL		30B
HTRF Human TNF alpha kit - Std		06AR
HTRF Human TNF alpha kit - 500 pts d2 antibody		09RA
HTRF Human TNF alpha kit - 500 pts Eu Cryptate antibody		09RA

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

hTNF alpha	Conc. (pg/ml)	DF %	DF Accepted Range Revision 2
Std 1	39	20%	13-33
Std 2	78	42%	34-61
Std 3	156	82%	70-119
Std 4	312	155%	120-216
Std 5	625	292%	220-359
Std 6	1250	530%	400-645
Std 7	2500	1068%	730-1100



Assay characteristics :

	Found value	Specifications Rev 02
Detection limit	6 pg/ml	< 39 pg/ml

HTRF Human TNF alpha kit - 500 pts - Lot n° : 09J

Accepted

Rejected

Quality control date : 2024-08-20	Supervisor's initials : FMV
-----------------------------------	-----------------------------

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.