

HTRF® Research Reagents

HTRF Human TNF alpha kit - 10K pts

Product Number : 62HTNFAPFH Lot Number : 11A

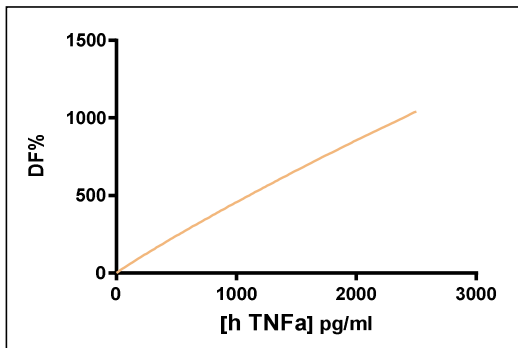
Expiration date : 2026-10-02 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 10 mL	21RA
HTRF Detect. Buf. 3 - 50 mL	29RA
HTRF Human TNF alpha kit - 10K pts d2 antibody	11AR
HTRF Human TNF alpha kit - 10K pts Eu Cryptate antibody	11AR
HTRF Human TNF alpha kit - Std	06AR
HTRF Human TNF alpha 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**

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	8 pg/ml	< 39 pg/ml

HTRF Human TNF alpha kit - 10K pts - Lot n° : 11A

Accepted

Rejected

Quality control date : 2023-12-28	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.