

HTRF® Research Reagents

HTRF Human TNF alpha kit - 10K pts

Product Number : 62HTNFAPEH Lot Number : 10B

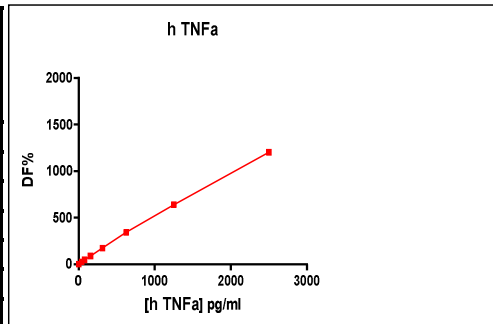
Expiration date : 2025-09-07 Document version : 1

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		10A
HTRF Human TNF alpha kit - 10K pts Eu Cryptate antibody		10
HTRF Human TNF alpha kit - Std		05RA
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	26%	13-33
Std 2	78	47%	34-61
Std 3	156	89%	70-119
Std 4	312	174%	120-216
Std 5	625	343%	220-359
Std 6	1250	641%	400-645
Std 7	2500	1203%	730-1100



*This batch has been validated despite Std 7 out of range

Assay characteristics :

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

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

Accepted

Rejected

Quality control date : 2023-10-03	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.