

HTRF® Research Reagents

HTRF Human TNF alpha kit - 500 pts

Product Number : 62HTNFAPEG Lot Number : 09G

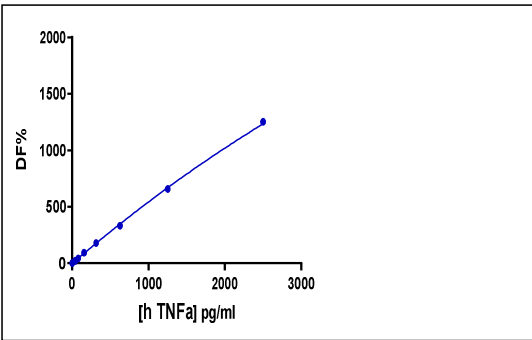
Expiration date : 2025-09-07 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 1.5 mL	30AR
HTRF Human TNF alpha kit - Std	05RA
HTRF Human TNF alpha kit - 500 pts d2 antibody	09RA
HTRF Human TNF alpha kit - 500 pts Eu Cryptate antibody	09RA
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	21%	13-33
Std 2	78	43%	34-61
Std 3	156	92%	70-119
Std 4	312	179%	120-216
Std 5	625	332%	220-359
Std 6	1250	657%	400-645
Std 7	2500	1252%	730-1100



This batch has been validated despite std 6 and 7 out of range

Assay characteristics :

	Found value	Specifications Rev 02
Detection limit	7 pg/ml	< 39 pg/ml

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

Accepted

Rejected

Quality control date : 2023-11-17	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.