

HTRF® Research Reagents

## HTRF Human TNF alpha kit - 96 pts

Product Number : 62HTNFAPET Lot Number : 09E

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

Reagents :	Lot n°
PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND	2203
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 0.5 mL	30AR
HTRF Human TNF alpha kit - Std	05RA
HTRF Human TNF alpha kit - 96 pts d2 antibody	09RA
HTRF Human TNF alpha kit - 96 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	30%	13-33
Std 2	78	55%	34-61
Std 3	156	102%	70-119
Std 4	312	200%	120-216
Std 5	625	359%	220-359
Std 6	1250	605%	400-645
Std 7	2500	915%	730-1100

**Assay characteristics :**

	Found value	Specifications Rev 02
Detection limit	5 pg/ml	≤ 39 pg/ml

HTRF Human TNF alpha kit - 96 pts - Lot n° : 09E

Accepted

Rejected

Quality control date : 2024-04-26	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.