

HTRF® Research Reagents

HTRF Human TNF alpha kit - 96 pts

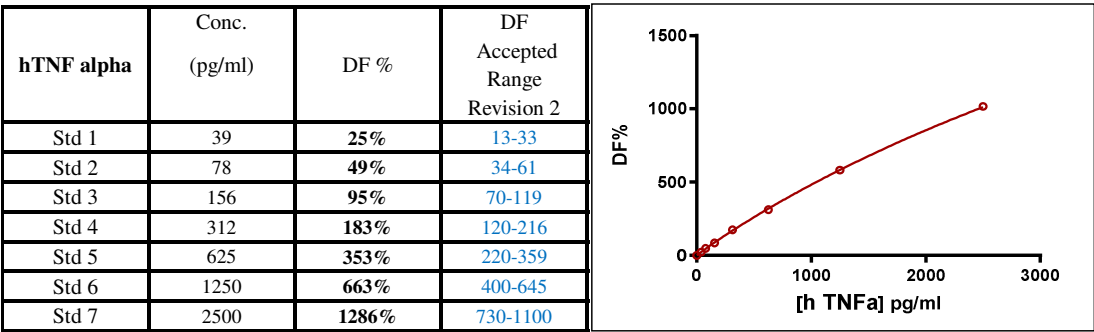
Product Number : 62HTNFAPET Lot Number : 09B

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	28RA
HTRF Human TNF alpha kit - Std	05A
HTRF Human TNF alpha kit - modop	8
HTRF Human TNF alpha kit - 96 pts d2 antibody	09A
HTRF Human TNF alpha kit - 96 pts Eu Cryptate antibody	09A

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**



Assay characteristics :

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

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

Accepted

Rejected

Quality control date : 2023-09-15	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 TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.