

HTRF® Research Reagents

HTRF Mouse TNF alpha kit - 500 pts

Product Number : 62MTNFAPEG Lot Number : 06C

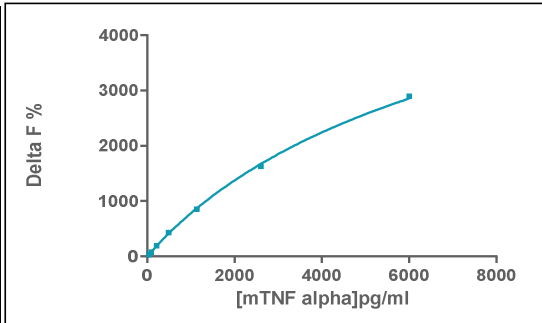
Expiration date : 2025-10-26 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 1.5 mL	30AR
HTRF Mouse TNF alpha kit - Std	06RA
HTRF Mouse TNF alpha kit - 500 pts d2 antibody	06RA
HTRF Mouse TNF alpha kit - 500 pts Eu Cryptate antibody	06RA
HTRF Mouse TNF alpha kit - modop	8

Results :

Incubation : 1h at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model : 4 Parameter Logistic (4PL 1/y2)* model

mTNFa	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	40.53	33%	25-46
Std 2	93.22	78%	61-102
Std 3	214.41	193%	143-228
Std 4	493.14	430%	212-493
Std 5	1134.22	855%	653-1029
Std 6	2608.7	1628%	1310-1980
Std 7	6000	2896%	2391-3271



Assay characteristics :

	Found value	Specifications Rev 01
Detection limite	4 pg/ml	≤ 40.5 pg/ml

HTRF Mouse TNF alpha kit - 500 pts - Lot n° : 06C

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.