

HTRF® Research Reagents

HTRF Hu. TNFa/TNFR1 Bind. kit - 500 pts

Product Number : 64BDTNFPEG Lot Number : 03D

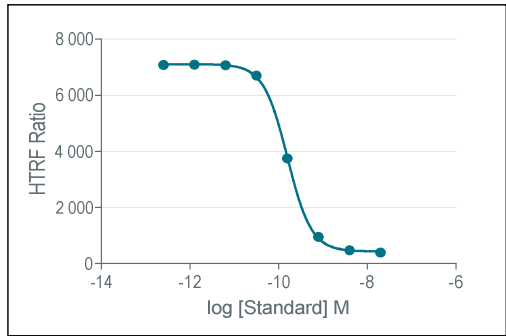
Expiration date : 2025-04-27 Document version : 1

Reagents :	Lot n°
HTRF PPI-Europium Detect. Buf. - 20 mL	05RA
HTRF Hu. TNFa/TNFR1 Bind. kit - 500 pts Tag2-TNFR1	03RA
HTRF Hu. TNFa/TNFR1 Bind. kit - Std	02RA
HTRF Hu. TNFa/TNFR1 Bind. kit - 500 pts Eu Cryptate antibody	03RB
HTRF Hu. TNFa/TNFR1 Bind. kit - 500 pts Tag1-Human TNF alpha	03RA
HTRF Hu. TNFa/TNFR1 Bind. kit - 500 pts XL665 reagent	03RA
HTRF Hu. TNFa/TNFR1 Bind. kit - modop	2

Results :

Incubation : Overnight at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model : log(inhibitor) vs. response -- Variable slope (four parameters)

	Conc. (nM)	RATIO
Std 0	0	7086
Std 1	0.128	7096
Std 2	0.32	7078
Std 3	0.64	6702
Std 4	1.6	3748
Std 5	8	951
Std 6	40	467
Std 7	200	392



Assay characteristics :

	Found value	Specifications Rev 01
ratio negatif	481	
S/B	18	
IC 50	0.16	0.15 - 0.27 nM

HTRF Hu. TNFa/TNFR1 Bind. kit - 500 pts - Lot n° : 03D Accepted Rejected

Quality control date : 2023-09-19 Supervisor's initials : FL

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.