

HTRF® Research Reagents

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

Product Number : 64BDTNFPEG Lot Number : 04C

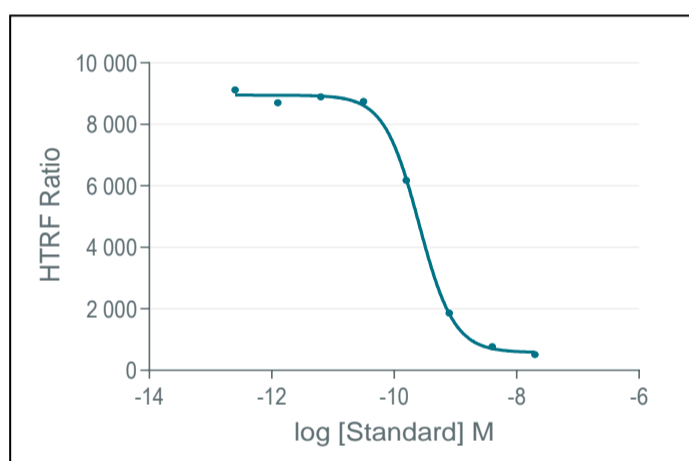
Expiration date : 2027-08-06 Document version : 1

Reagents :	Lot n°
HTRF PPI-Europium Detect. Buf. - 20 mL	08A
HTRF Hu. TNFa/TNFR1 Bind. kit - 500 pts Tag2-TNFR1	04B
HTRF Hu. TNFa/TNFR1 Bind. kit - Std	05A
HTRF Hu. TNFa/TNFR1 Bind. kit - 500 pts Eu Cryptate antibody	04A
HTRF Hu. TNFa/TNFR1 Bind. kit - 500 pts Tag1-Human TNF alpha	04B
HTRF Hu. TNFa/TNFR1 Bind. kit - 500 pts XL665 reagent	04A

Results :

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

	Conc. (nM)	RATIO
Std 0	0	9120
Std 1	0,128	8700
Std 2	0,32	8894
Std 3	0,64	8748
Std 4	1,6	6175
Std 5	8	1866
Std 6	40	773
Std 7	200	516



Assay characteristics :

	Found value	Specifications Rev 01
Negative	700	
S/B	18	
IC 50	0,25	0.15-0.27 nM

HTRF Hu. TNFa/TNFR1 Bind. kit - 500 pts - Lot n° : 04C

Accepted

Rejected

Quality control date :	2025-01-03	Supervisor's initials :	MP
------------------------	------------	-------------------------	----

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.