

HTRF® Research Reagents

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

Product Number : 64BDTNFPEG Lot Number : 05A

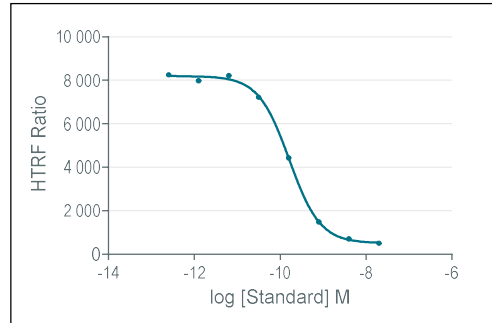
Expiration date : 2027-09-09 Document version : 1

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

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	8252
Std 1	0,128	7981
Std 2	0,32	8225
Std 3	0,64	7223
Std 4	1,6	4430
Std 5	8	1485
Std 6	40	700
Std 7	200	502



Assay characteristics :

	Found value	Specifications Rev 01
Negative	706	
S/B	16	
IC 50	0,163	0.15-0.27 nM

HTRF Hu. TNFa/TNFR1 Bind. kit - 500 pts - Lot n° : 05A

Accepted

Rejected

Quality control date : 2024-09-18	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.