

HTRF® Research Reagents

HTRF IL17A/IL17RA Bind. kit - 500 pts

Product Number : 64BDIL17PEG Lot Number : 06G

Expiration date : 2025-10-18 Document version : 1

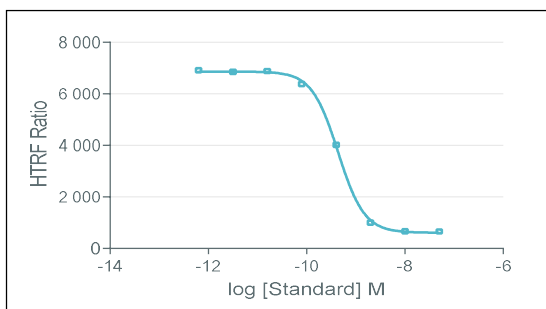
Reagents :	Lot n°
HTRF PPI-Europium Detect. Buf. - 20 mL	07RA
HTRF IL17A/IL17RA Bind. kit - 500 pts Tag2-Human IL17RA	09AR
HTRF IL17A/IL17RA Bind. kit - Std	04RA
HTRF IL17A/IL17RA Bind. kit - 500 pts Eu Cryptate reagent	06RA
HTRF IL17A/IL17RA Bind. kit - 500 pts Tag1-Human IL17A	09AR
HTRF IL17A/IL17RA Bind. kit - 500 pts XL665 antibody	06B

Results :

Incubation : Overnight at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model : nonlinear regression, sigmoidal dose response (variable slope)

Assay characteristics :

hIL17A/IL17RA	Conc. (pMi)	RATIO
Std 0	0	6917
Std 1	3,2000	6855
Std 2	16,000	6880
Std 3	80,000	6383
Std 4	400,0	4018
Std 5	2000	1011
Std 6	10000	672
Std 7	50000	660



Assay characteristics :

	Found value	Specifications Rev 01
negative control (ratio)	433	350 - 531
IC 50	0,44	0.3 - 0.6
S/B	10	8 - 14

HTRF IL17A/IL17RA Bind. kit - 500 pts - Lot n° : 06G

Accepted

Rejected

Quality control date : 2024-06-18	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.