

HTRF® Research Reagents

HTRF IL6/IL6R Bind. kit - 500 pts

Product Number : 64BDIL6PEG Lot Number : 01C

Expiration date : 2025-08-06 Document version : 1

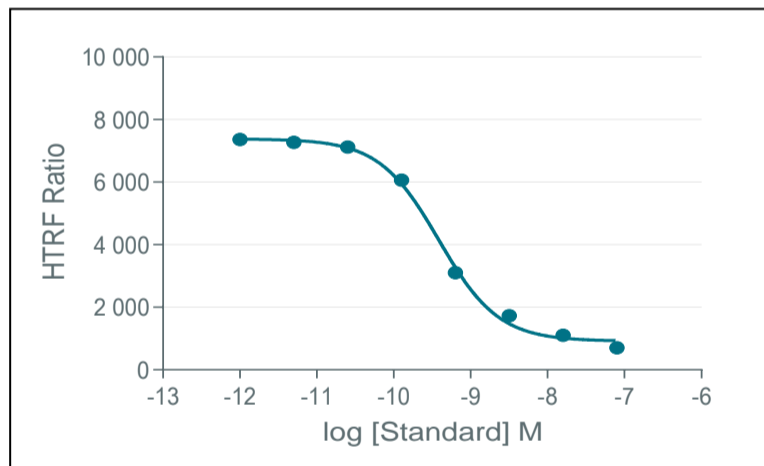
Reagents :	Lot n°
HTRF PPI-Europium Detect. Buf. - 20 mL	07RA
HTRF IL6/IL6R Bind. kit - 500 pts Tag2-Human IL6R	01RB
HTRF IL6/IL6R Bind. kit - Std	01RA
HTRF IL6/IL6R Bind. kit - 500 pts Eu Cryptate antibody	01RB
HTRF IL6/IL6R Bind. kit - 500 pts Tag1-Human IL6	01RB
HTRF IL6/IL6R Bind. kit - 500 pts XL665 antibody	01RB

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	7360
Std 1	3,2000	7273
Std 2	16,000	7120
Std 3	80,000	6062
Std 4	400,0	3100
Std 5	2000	1727
Std 6	10000	1101
Std 7	50000	702



Assay characteristics :

	Found value	Specifications
negative control (ratio)	497	488 - 564
IC 50	0,38	0.14 -0.62
S/B	10	

HTRF IL6/IL6R Bind. kit - 500 pts - Lot n° : 01C

Accepted

Rejected

Quality control date : 2024-02-09	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.