

HTRF® Research Reagents

HTRF Human IL17 kit - 500 pts

Product Number : 62HIL17PEG Lot Number : 04E

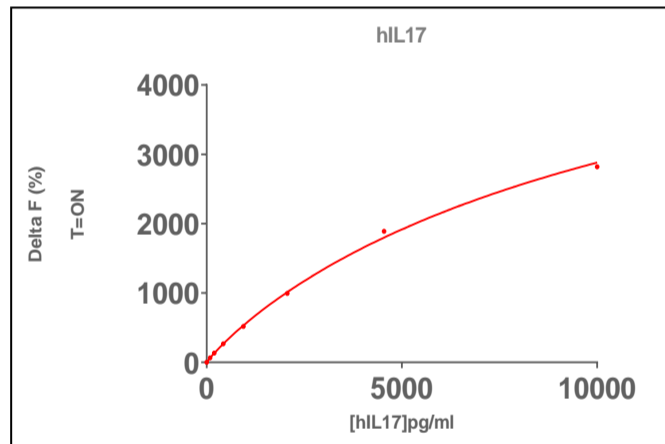
Expiration date : 2027-04-24 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21B
HTRF Detect. Buf. 3 - 1.5 mL	30B
HTRF Human IL17 kit - Std	04A
HTRF Human IL17 kit - 500 pts d2 antibody	04B
HTRF Human IL17 kit - 500 pts Eu Cryptate antibody	04C

Results :

Incubation : Overnight at RT  
 Plate : 384 wells small volume white plate 20 µL  
 Reader : Pherastar Plus  
 Fitting model : 4 Parameter Logistic (4PL 1/y2)\* model

IL17	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	88	64%	48-66
Std 2	194	134%	99-134
Std 3	427	268%	185-275
Std 4	939	518%	366-612
Std 5	2066	993%	739-1275
Std 6	4545	1890%	1556-2551
Std 7	10000	2820%	2676-3186



Assay characteristics :

	Found value	Specifications Rev 02
Detection limit	5,5 pg/ml	≤ 8 pg/ml

HTRF Human IL17 kit - 500 pts - Lot n° : 04E

Accepted

Rejected

Quality control date : 2024-07-01	Supervisor's initials : PL
-----------------------------------	----------------------------

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.