

HTRF® Research Reagents

HTRF Human IL17 kit - 96 pts

Product Number : 62HIL17PET Lot Number : 01P

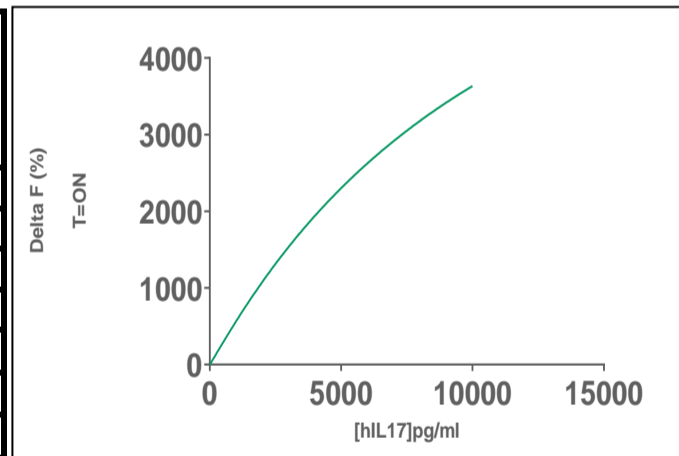
Expiration date : 2025-11-05 Document version : 1

Reagents :	Lot n°
PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND	2203
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 0.5 mL	28RA
HTRF Human IL17 kit - Std	03RA
HTRF Human IL17 kit - 96 pts d2 antibody	01RA
HTRF Human IL17 kit - 96 pts Eu Cryptate antibody	01RA

Results :

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

IL17	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	88	52%	48-66
Std 2	194	105%	99-134
Std 3	427	227%	185-275
Std 4	939	505%	366-612
Std 5	2066	1123%	739-1275
Std 6	4545	2420%	1556-2551
Std 7	10000	3383%	2676-3186



\* This batch has been validated despite Std 7 out of range

Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	6,8 pg/ml	≤ 8pg/ml

HTRF Human IL17 kit - 96 pts - Lot n° : 01P

Accepted

Rejected

Quality control date : 2024-04-15	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.