

HTRF® Research Reagents

HTRF Human IL17 kit - 10K pts

Product Number : 62HIL17PEH Lot Number : 03E

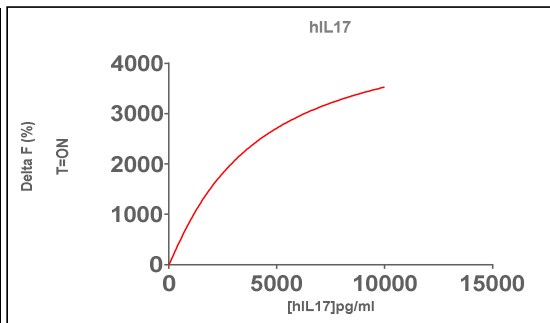
Expiration date : 2024-07-21 Document version : 1

Reagents :		Lot n°
HTRF Diluent 5 (5X) - 10 mL		21RA
HTRF Detect. Buf. 3 - 50 mL		29RA
HTRF Human IL17 kit - 10K pts d2 antibody		03RA
HTRF Human IL17 kit - 10K pts Eu Cryptate antibody		03RA
HTRF Human IL17 kit - Std		03RA
HTRF Human IL17 kit - modop		7

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	79%	39-52
Std 2	194	176%	95-110
Std 3	427	378%	199-245
Std 4	939	785%	444-522
Std 5	2066	1665%	924-1194
Std 6	4545	2832%	1940-2485
Std 7	10000	3274%	2938-3363



Assay characteristics :

	Found value	Specifications Rev 02
Detection limite	4.5 pg/ml	≤ 8 pg/ml

HTRF Human IL17 kit - 10K pts - Lot n° : 03E

Accepted

Rejected

Quality control date :	2023-11-09	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 TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.