

HTRF® Research Reagents

HTRF Mouse IL17 kit - 500 pts

Product Number : 62MIL17PEG Lot Number : 04J

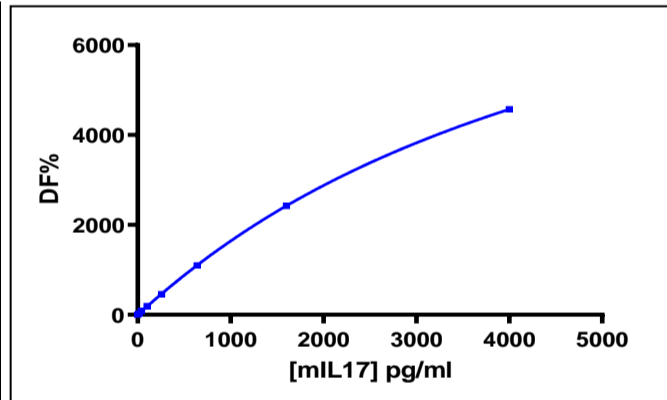
Expiration date : 2025-07-06 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21B
HTRF Detect. Buf. 3 - 1.5 mL	30B
HTRF Mouse IL17 kit - Std	02RA
HTRF Mouse IL17 kit - 500 pts d2 antibody	04RA
HTRF Mouse IL17 kit - 500 pts Eu Cryptate antibody	04RA

Results :

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

mIL17	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	16	33%	25-59
Std 2	41	80%	68-139
Std 3	102	191%	170-314
Std 4	256	459%	427-739
Std 5	640	1099%	1007-1653
Std 6	1600	2423%	2200-3246
Std 7	4000	4571%	3507-4615



Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	2 pg/ml	≤ 16 pg/ml

HTRF Mouse IL17 kit - 500 pts - Lot n° : 04J

Accepted

Rejected

Quality control date : 2024-12-05	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.