

HTRF® Research Reagents

# HTRF Human CCL17 Kit - 500 pts

Product Number : 62HCCL17PEG Lot Number : 01A

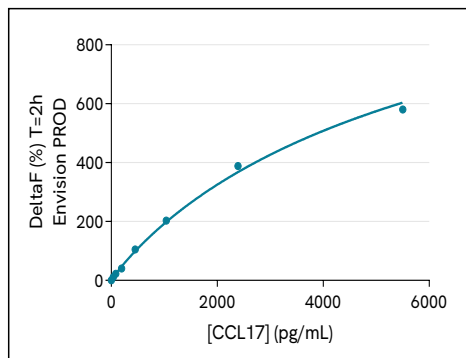
Expiration date : 2027-12-15 Document version :

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21B
HTRF Detect. Buf. 3 - 1.5 mL	30B
HTRF Human CCL17 Kit - Std	01A
HTRF Human CCL17 Kit - 500 pts d2 antibody	01A
HTRF Human CCL17 Kit - 500 pts Eu Cryptate antibody	01A

**Results :**

Incubation : 2h at RT  
 Plate : 384 wells small volume white plate 20 µL  
 Reader : EnVision  
 Fitting model : 4 Parameter Logistic (4PL)

hCCL17	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	37,2	14%	11- 19
Std 2	85,5	32%	23-35
Std 3	197	71%	46-73
Std 4	452	150%	112 - 168
Std 5	1040	296%	214-321
Std 6	2391	553%	398-598
Std 7	5500	595%	469 - 703



**Assay characteristics :**

	Found value	Specifications Rev 01
detection limite	9,7	< 37 pg/ml

HTRF Human CCL17 Kit - 500 pts - Lot n° : 01A

Accepted

Rejected

Quality control date : 2025-04-07	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.