

HTRF® Research Reagents

HTRF Human CXCL1 kit - 500 pts

Product Number : 62HCXC1PEG Lot Number : 06E

Expiration date : 2026/03/29 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 1.5 mL	29A
HTRF Human CXCL1 kit - Std	05AR
HTRF Human CXCL1 kit - 500 pts d2 antibody	06A
HTRF Human CXCL1 kit - 500 pts Eu Cryptate antibody	06A
HTRF Human CXCL1 kit - modop	7

Results :

Incubation : 2h at 4°C
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model : 4 Parameter Logistic (4PL 1/y2)* model

CXCL 1	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	118	52%	45-75
Std 2	212	107%	80-140
Std 3	381	179%	150-230
Std 4	686	291%	235-350
Std 5	1235	419%	360-470
Std 6	2222	574%	480-600
Std 7	4000	705%	606-780

Assay characteristics :

	Found value	Specifications Rev 02
Detection limit	4 pg/ml	≤ 10 pg/ml
	/	/
	/	/
	/	/

HTRF Human CXCL1 kit - 500 pts - Lot n° : 06E

Accepted

Rejected

Quality control date : 2023-09-26	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.