

HTRF® Research Reagents

HTRF Human CXCL1 kit - 500 pts

Product Number : 62HCXC1PEG Lot Number : 06C

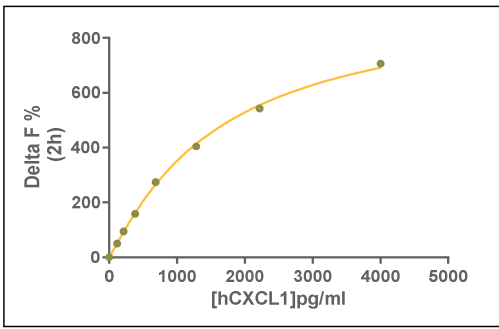
Expiration date : 2025-05-01 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 1.5 mL	28RA
HTRF Human CXCL1 kit - Std	04RA
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	50%	45-75
Std 2	212	94%	80-140
Std 3	381	159%	150-230
Std 4	686	274%	235-350
Std 5	1235	404%	360-470
Std 6	2222	543%	480-600
Std 7	4000	706%	606-780



Assay characteristics :

	Found value	Specifications Rev 02
Detection limit	9.4 pg/ml	≤ 10 pg/ml
	/	/
	/	/
	/	/

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

Accepted

Rejected

Quality control date : 2023-09-01 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.