

HTRF® Research Reagents

HTRF Human CXCL1 kit - 10K pts

Product Number : 62HCXC1PEH Lot Number : 08A

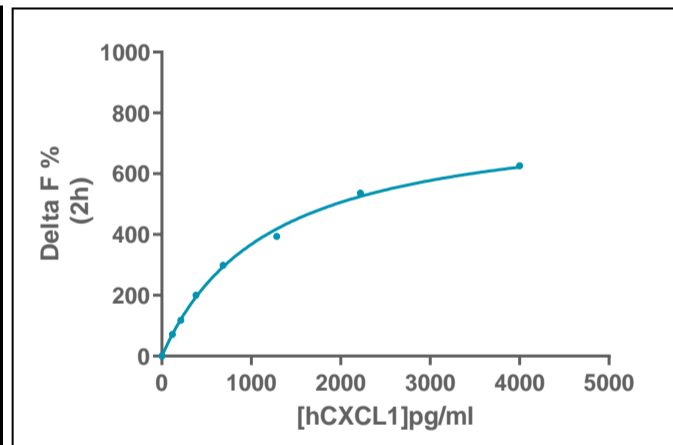
Expiration date : 2026-05-23 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 10 mL	21RA
HTRF Detect. Buf. 3 - 50 mL	29RA
HTRF Human CXCL1 kit - 10K pts d2 antibody	08A
HTRF Human CXCL1 kit - 10K pts Eu Cryptate antibody	08A
HTRF Human CXCL1 kit - Std	05AR

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 Revision n°2
Std 1	118	71%	45-75
Std 2	212	118%	80-140
Std 3	381	200%	150-230
Std 4	686	299%	235-350
Std 5	1235	394%	360-470
Std 6	2222	536%	480-600
Std 7	4000	626%	606-780



Assay characteristics :

	Found value	Specifications Rev 02
Detection limit	8.5 pg/ml	≤ 10 pg/ml

HTRF Human CXCL1 kit - 10K pts - Lot n° : 08A

Accepted

Rejected

Quality control date : 2024-07-03	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.