

HTRF® Research Reagents

HTRF Human CXCL10 kit - 10K pts

Product Number : 62HCX10PEH Lot Number : 10A

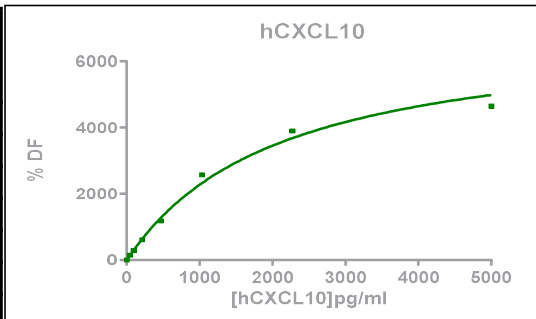
Expiration date : 2024-09-01 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 10 mL	21RA
HTRF Detect. Buf. 3 - 50 mL	29RA
HTRF Human CXCL10 kit - 10K pts d2 antibody	10A
HTRF Human CXCL10 kit - 10K pts Eu Cryptate antibody	10A
HTRF Human CXCL10 kit - Std	05RA

Results :

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

hCXCL10	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	44,1	147%	62 --222
Std 2	97	287%	225-446
Std 3	213,4	608%	450-880
Std 4	496,6	1178%	900-1481
Std 5	1033,1	2574%	1599-2649
Std 6	2272,7	3895%	2828-4345
Std 7	5000	4649%	3447-5287



Assay characteristics :

	Found value	Specifications Rev 03
Detection limit	1.3 pg/ml	≤ 44.1 pg/ml

HTRF Human CXCL10 kit - 10K pts - Lot n° : 10A

Accepted

Rejected

Quality control date : 2024-02-19	Supervisor's initials : MP
-----------------------------------	----------------------------

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.