

HTRF® Research Reagents

HTRF Human CCL22 Kit - 500 pts

Product Number : 62HCCL22PEG Lot Number : 01C

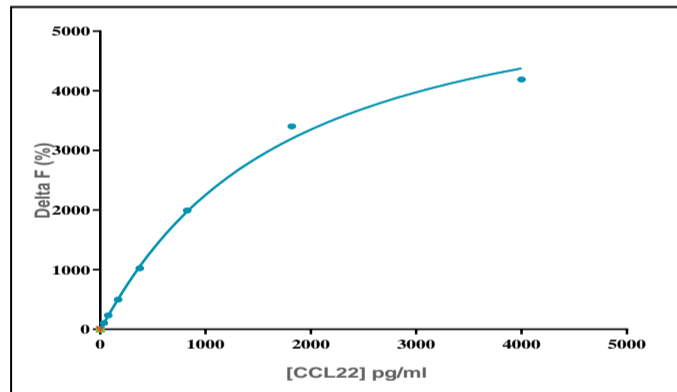
Expiration date : 2025-07-19 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21B
HTRF Detect. Buf. 3 - 1.5 mL	30B
HTRF Human CCL22 Kit - Std	01RA
HTRF Human CCL22 Kit - 500 pts d2 antibody	01RA
HTRF Human CCL22 Kit - 500 pts Eu Cryptate antibody	01RA

Results :

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

hCCL22	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	35	105%	84 -- 124
Std 2	78	233%	160 --- 306
Std 3	171	497%	348 --- 646
Std 4	376	1023%	800 --- 1247
Std 5	826	1993%	1594 -- 2392
Std 6	1818	3403%	2722 -- 4084
Std 7	4000	4190%	3352 -- 5028



Assay characteristics :

	Found value	Specifications Rev 02
Detection limit	1 pg/ml	< 35 pg/ml

HTRF Human CCL22 Kit - 500 pts - Lot n° : 01C

Accepted

Rejected

Quality control date : 2024-07-30	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.