

HTRF® Research Reagents

HTRF Human CCL4 kit - 10K pts

Product Number : 62HCCL4PEH Lot Number : 05D

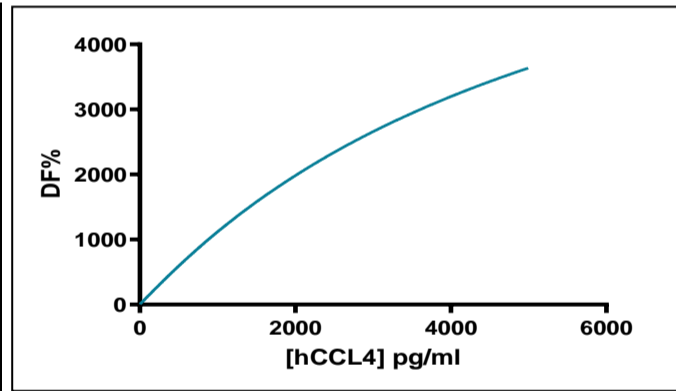
Expiration date : 2025-07-17 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 10 mL	21RA
HTRF Detect. Buf. 3 - 50 mL	29RA
HTRF Human CCL4 kit - 10K pts d2 antibody	05RA
HTRF Human CCL4 kit - 10K pts Eu Cryptate antibody	05RA
HTRF Human CCL4 kit - Std	04RA

Results :

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

hCCL2	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	33,8	46%	33-52
Std 2	77,7	104%	62-118
Std 3	178,7	223%	140-261
Std 4	411	487%	327-599
Std 5	945,2	1064%	732-1264
Std 6	2173,9	2113%	1555-2406
Std 7	5000	3636%	2778-3810



Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	2.8 pg/ml	≤ 33.8 pg/ml

HTRF Human CCL4 kit - 10K pts - Lot n° : 05D

Accepted

Rejected

Quality control date : 2024-04-22	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.