

HTRF® Research Reagents

HTRF Human CCL2 kit - 500 pts

Product Number : 62HCCL2PEG Lot Number : 06H

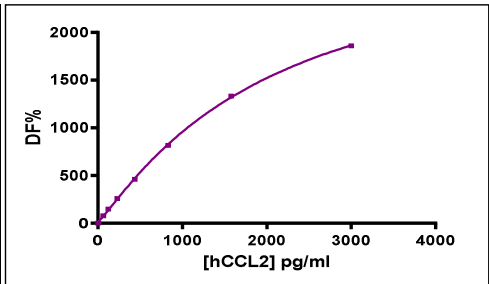
Expiration date : 2025-05-27 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 1.5 mL	29A
HTRF Human CCL2 kit - Std	05RA
HTRF Human CCL2 kit - 500 pts d2 antibody	06RA
HTRF Human CCL2 kit - 500 pts Eu Cryptate antibody	06RA
HTRF Human CCL2 kit - modop	7

Results :

Incubation : Overnight 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	63.77	78%	50-106
Std 2	121.16	148%	116-199
Std 3	230.2	258%	220-355
Std 4	437.38	461%	361-616
Std 5	831.02	816%	653-988
Std 6	1578.95	1334%	1074-1410
Std 7	3000	1860%	1440-1910



Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	3.3 pg/ml	≤ 8 pg/ml
	/	/
	/	/
	/	/

HTRF Human CCL2 kit - 500 pts - Lot n° : 06H

Accepted

Rejected

Quality control date : 2023-08-16	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.