

HTRF® Research Reagents

HTRF Human CCL3 kit - 500 pts

Product Number : 62HCCL3PEG Lot Number : 02E

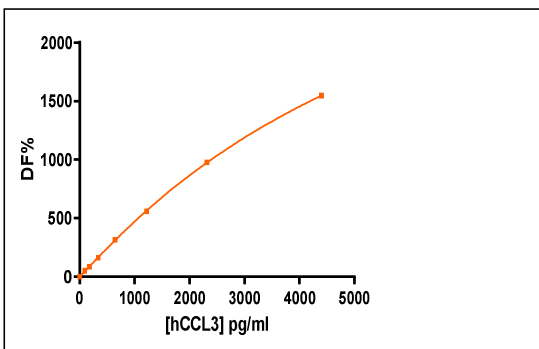
Expiration date : 2024-07-28 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 1.5 mL	30AR
HTRF Human CCL3 kit - Std	02RB
HTRF Human CCL3 kit - 500 pts d2 antibody	02RA
HTRF Human CCL3 kit - 500 pts Eu Cryptate antibody	02RA

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

Hccl3	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	93.5	53%	22-57
Std 2	177.7	85%	57-86
Std 3	337.6	163%	119-178
Std 4	641.5	316%	220-330
Std 5	1218.8	559%	445-668
Std 6	2315.8	978%	874-1311
Std 7	4400.0	1547%	1513-2269



Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	7 pg/ml	≤ 93.5 pg/ml

HTRF Human CCL3 kit - 500 pts - Lot n° : 02E

Accepted

Rejected

Quality control date : 2023-11-29	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.