

HTRF® Research Reagents

HTRF Mouse CCL2 kit - 500 pts

Product Number : 62MCCL2PEG Lot Number : 03G

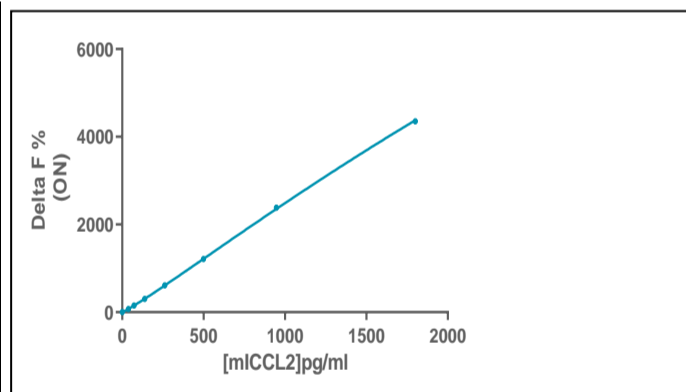
Expiration date : 2025-11-27 Document version : 1

Reagents :		Lot n°
HTRF Diluent 5 (5X) - 2 mL		21RA
HTRF Detect. Buf. 3 - 1.5 mL		30AR
HTRF Mouse CCL2 kit - Std		02RA
HTRF Mouse CCL2 kit - 500 pts d2 antibody		03B
HTRF Mouse CCL2 kit - 500 pts Eu Cryptate antibody		03B

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

mCCL2	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	38,3	74%	53-69
Std 2	72,7	152%	113-149
Std 3	138,1	300%	206-341
Std 4	262,4	612%	397-746
Std 5	498,6	1212%	773-1567
Std 6	947,4	2381%	1554-3307
Std 7	1800	4353%	3276-5875



Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	9 pg/ml	< 38 pg/ml

HTRF Mouse CCL2 kit - 500 pts - Lot n° : 03G

Accepted

Rejected

Quality control date : 2024-07-03	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.