

HTRF® Research Reagents

HTRF Human IL12 kit - 500 pts

Product Number : 62HIL12PEG Lot Number : 05F

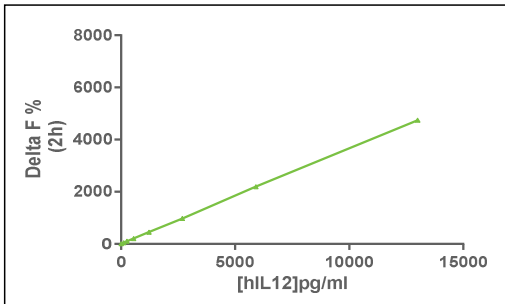
Expiration date : 2025-06-27 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 1.5 mL	29A
HTRF Human IL12 kit - Std	07RA
HTRF Human IL12 kit - 500 pts d2 antibody	05RA
HTRF Human IL12 kit - 500 pts Eu Cryptate antibody	05RA
HTRF Human IL12 kit - modop	7

Results :

Incubation : 2h at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model : 4 Parameter Logistic (4PL 1/y2)* model

hIL12	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	115	48%	24-69
Std 2	252	106%	60-153
Std 3	555	211%	150-296
Std 4	1221	452%	341-651
Std 5	2686	978%	775-1440
Std 6	5909	2196%	1830-3219
Std 7	13000	4742%	4194-6643



Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	10 pg/ml	< 115 pg/ml
	/	/
	/	/
	/	/

HTRF Human IL12 kit - 500 pts - Lot n° : 05F

Accepted

Rejected

Quality control date : 2023-10-03	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.