

HTRF® Research Reagents

HTRF Human IL12 kit - 10K pts

Product Number : 62HIL12PEH Lot Number : 07B

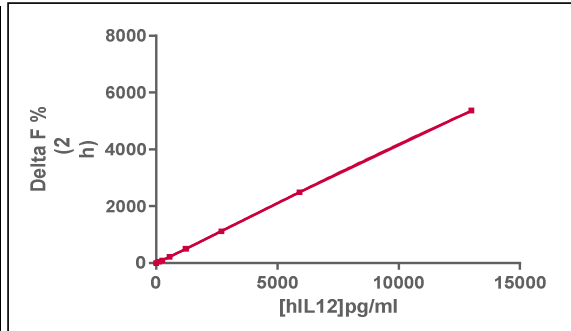
Expiration date : 2025-11-07 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 10 mL	21RA
HTRF Detect. Buf. 3 - 50 mL	29RA
HTRF Human IL12 kit - 10K pts d2 antibody	07A
HTRF Human IL12 kit - 10K pts Eu Cryptate antibody	07A
HTRF Human IL12 kit - Std	07RA

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	42%	24-69
Std 2	252	95%	60-153
Std 3	555	218%	150-296
Std 4	1221	503%	341-651
Std 5	2686	1117%	775-1440
Std 6	5909	2487%	1830-3219
Std 7	13000	5368%	4194-6643



Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	11 pg/ml	< 115 pg/ml

HTRF Human IL12 kit - 10K pts - Lot n° : 07B

Accepted

Rejected

Quality control date : 2023-11-30	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.