

HTRF® Research Reagents

HTRF Human IL22 kit - 500 pts

Product Number : 62HIL22PEG Lot Number : 03D

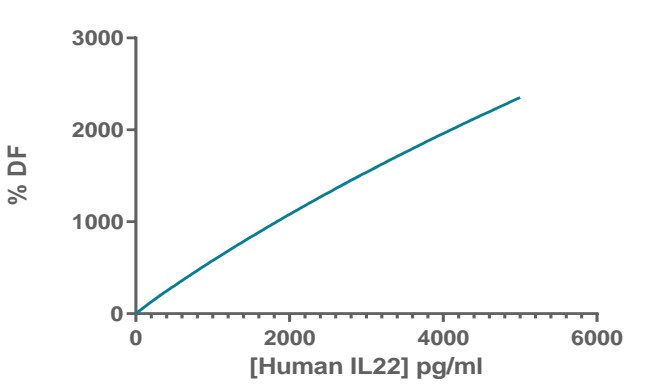
Expiration date : 2025-12-19 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21B
HTRF Detect. Buf. 3 - 1.5 mL	30B
HTRF Human IL22 kit - Std	02RA
HTRF Human IL22 kit - 500 pts d2 antibody	03RA
HTRF Human IL22 kit - 500 pts Eu Cryptate antibody	03RA

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

hIL22	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	34	24%	17 -- 26
Std 2	78	51%	46 -- 69
Std 3	179	114%	92 -- 139
Std 4	411	241%	218 -- 327
Std 5	945	573%	472 -- 708
Std 6	2174	1165%	926 -- 1389
Std 7	5000	2335%	1577 -- 2365



Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	6 pg/ml	< 34 pg/ml

HTRF Human IL22 kit - 500 pts - Lot n° : 03D

Accepted

Rejected

Quality control date : 2024-07-02	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.