

HTRF® Research Reagents

HTRF Human IL4 kit - 500 pts

Product Number : 62HIL04PEG Lot Number : 02S

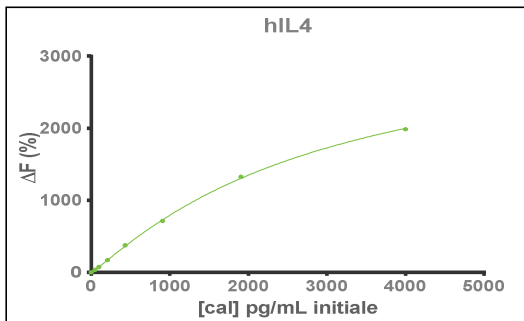
Expiration date : 2024-12-06 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	19G
HTRF Detect. Buf. 3 - 1.5 mL	28A
HTRF Human IL4 kit - Std	04RA
HTRF Human IL4 kit - 500 pts d2 antibody	02RC
HTRF Human IL4 kit - 500 pts Eu Cryptate antibody	02RC
HTRF Human IL4 kit - modop	7

Results :

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

IL4	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	47	35%	35-63
Std 2	98	74%	59-132
Std 3	206	169%	128-257
Std 4	432	375%	249-520
Std 5	907	713%	489-959
Std 6	1905	1324%	880-1678
Std 7	4000	1984%	1583-2106



Assay characteristics :

	Found value	Specifications Rev 01
Detection limite	6 pg/ml	≤ 47 pg/ml

HTRF Human IL4 kit - 500 pts - Lot n° : 02S

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.