

HTRF® Research Reagents

HTRF Human IL2 kit - 96 pts

Product Number : 62HIL02PET

Lot Number : 01ZF

Expiration date : 2026-03-20

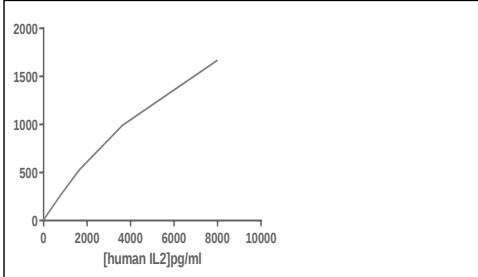
Document version : 1

Reagents :		Lot n°
PLAQUE BLANCHE 96 PUIITS SV		2450
PLAQUE BLANCHE 96 PUIITS SV		2450
HTRF Diluent 5 (5X) - 2 mL		22A
HTRF Detect. Buf. 3 - 0.5 mL		30AR
HTRF Human IL2 kit - Std		07RA
HTRF Human IL2 kit - 96 pts d2 antibody		01RA
HTRF Human IL2 kit - 96 pts Eu Cryptate antibody		01RA

Results :

Incubation : **3h at RT**
Plate : **384 wells small volume white plate 20 µL**
Reader : **Pherastar**
Fitting model : **4 Parameter Logistic (4PL 1/y2)* model**

IL2	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	70,56	27%	10-45
Std 2	155,23	58%	35-89
Std 3	341,51	121%	100-172
Std 4	751,3	255%	182-363
Std 5	1652,9	530%	389-695
Std 6	3636,4	993%	742-1171
Std 7	8000	1670%	1172-1670



Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	11 pg/ml	≤ 20 pg/ml

HTRF Human IL2 kit - 96 pts - Lot n° : 01ZF

☒ Accepted☐ Rejected

Quality control date : 2025-04-14

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 TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.