

## HTRF® Research Reagents

## HTRF Human IL2 kit - 10K pts

Product Number : 62HIL02PEH

Lot Number : 07F

Expiration date : 2024-11-25

Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 10 mL	21A
HTRF Detect. Buf. 3 - 50 mL	29A
HTRF Human IL2 kit - 10K pts d2 antibody	07RA
HTRF Human IL2 kit - 10K pts Eu Cryptate antibody	07RA
HTRF Human IL2 kit - Std	07A
HTRF Human IL2 kit - modop	7

## Results :

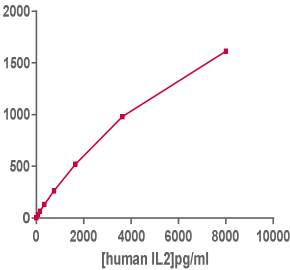
Incubation : 3h at RT

Plate : 384 wells small volume white plate 20 µL

Reader : Pherastar Plus

Fitting model : 4 Parameter Logistic (4PL 1/y2)\* model

IL2	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	70.56	31%	10-45
Std 2	155.23	65%	35-89
Std 3	341.51	130%	100-172
Std 4	751.3	261%	182-363
Std 5	1652.9	521%	389-695
Std 6	3636.4	980%	742-1171
Std 7	8000	1613%	1172-1670



## Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	9 pg/ml	< 70.6 pg/ml
	/	/
	/	/
	/	/

HTRF Human IL2 kit - 10K pts - Lot n° : 07F

☒ Accepted☐ Rejected

Quality control date : 2023-08-11

Supervisor's initials : MP

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.