

HTRF® Research Reagents

HTRF Human IL10 kit - 500 pts

Product Number :62HIL10PEG

Lot Number :06J

Expiration date :2026-11-17

Document version :1

Reagents :		Lot n°
HTRF Diluent 5 (5X) - 2 mL		21B
HTRF Detect. Buf. 3 - 1.5 mL		30B
HTRF Human IL10 kit - Std		06RA
HTRF Human IL10 kit - 500 pts d2 antibody		06RA
HTRF Human IL10 kit - 500 pts Eu Cryptate antibody		06RA

Results :

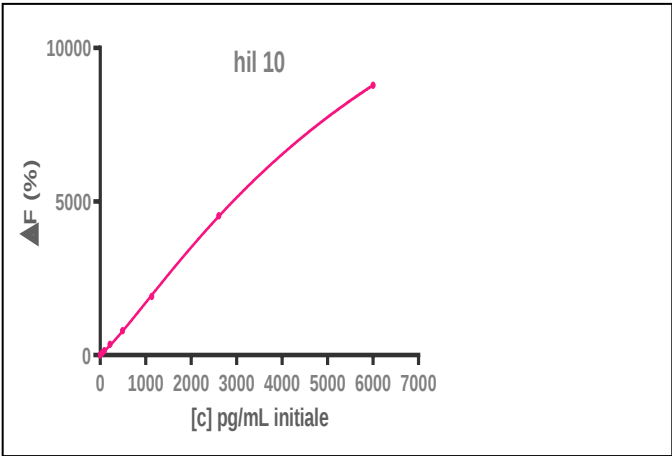
Incubation :Overnight at RT

Plate :384 wells small volume white plate 20 µL

Reader :Pherastar Plus

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

IL10	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	40	68%	27-77
Std 2	93	146%	73-185
Std 3	214	351%	203-433
Std 4	493	797%	608-1028
Std 5	1134	1913%	1642-2614
Std 6	2608	4541%	4395-7131
Std 7	6000	8786%	8744-10139



Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	2pg/ml	≤ 40 pg/ml

HTRF Human IL10 kit - 500 pts - Lot n° : 06J

☒ Accepted

☐ Rejected

Quality control date :2024-08-20	Supervisor's initials :FMV
----------------------------------	----------------------------

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.