

HTRF® Research Reagents

HTRF Human IL10 kit - 500 pts

Product Number :62HIL10PEG

Lot Number :06L

Expiration date : 2027-03-17

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		07A
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	73%	27-77	
Std 2	93	178%	73-185	
Std 3	214	415%	203-433	
Std 4	493	968%	608-1028	
Std 5	1134	2510%	1642-2614	
Std 6	2608	6533%	4395-7131	
Std 7	6000	9900%	8744-10139	

Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	2,2 pg/ml	≤ 40 pg/ml
HTRF Human IL10 kit - 500 pts - Lot n° : 06L		
☒ Accepted ☐ Rejected		
Quality control date :2024-12-06	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.