

HTRF® Research Reagents

HTRF Human IL11 kit - 500 pts

Product Number : 62HIL11PEG Lot Number : 01F

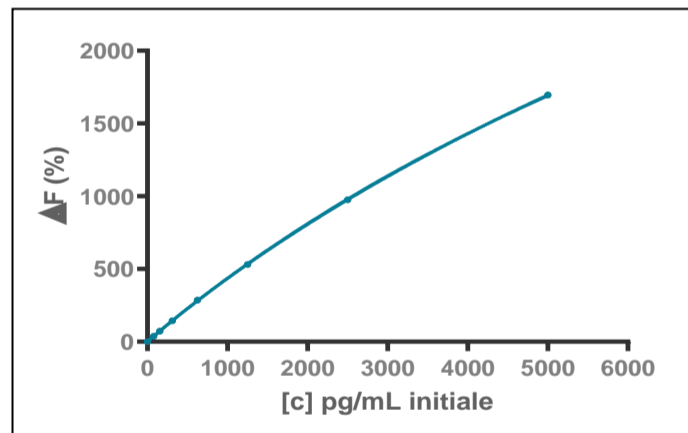
Expiration date : 2026-10-27 Document version : 1

Reagents :		Lot n°
HTRF Diluent 5 (5X) - 2 mL		21B
HTRF Detect. Buf. 3 - 1.5 mL		30B
HTRF Human IL11 kit - Std		02RA
HTRF Human IL11 kit - 500 pts d2 antibody		01RA
HTRF Human IL11 kit - 500 pts Eu Cryptate antibody		01RA

Results :

Incubation : Overnight at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model : 4 Parameter Logistic (4PL)

h IL11	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	78,1	CV	12-92
Std 2	156,3	0%	47-149
Std 3	312,5	0%	101-266
Std 4	625,0	0%	218-441
Std 5	1250,0	0%	456-801
Std 6	2500,0	0%	879-1499
Std 7	5000,0	0%	1674-2920



Assay characteristics :

	Found value	Specifications
Detection limite	8 pg/ml	< 78 pg/ml

HTRF Human IL11 kit - 500 pts - Lot n° : 01F

Accepted

Rejected

Quality control date : 2024-09-25	Supervisor's initials : FL
-----------------------------------	----------------------------

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.