

HTRF® Research Reagents

HTRF Human IL11 kit - 500 pts

Product Number : 62HIL11PEG **Lot Number :** 01E

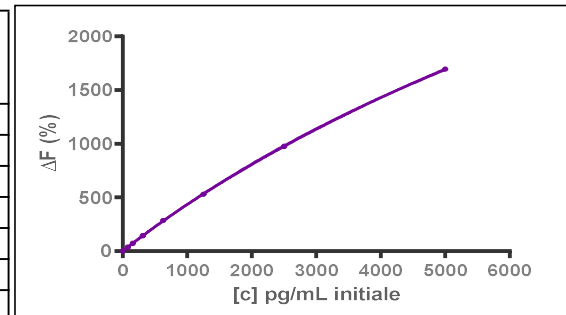
Expiration date : 2026-10-27 **Document version :** 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 1.5 mL	30AR
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)

M IFN gamma	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	78.1	REV n°1	12-92
Std 2	156.3	0%	47-149
Std 3	312.5	12-92	101-266
Std 4	625.0	47-149	218-441
Std 5	1250.0	101-266	456-801
Std 6	2500.0	218-441	879-1499
Std 7	5000.0	456-801	1674-2920


Assay characteristics :

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

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

 Rejected

Quality control date : 2022-04-14 **Supervisor's initials :** AC

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.