

HTRF® Research Reagents

HTRF Mouse IFN g kit - 500 pts

Product Number :

62MIFNGPEG

Lot Number :

05G

Expiration date :

2026-01-26

Document version :

1

Reagents :		Lot n°
HTRF Diluent 5 (5X) - 2 mL		21B
HTRF Detect. Buf. 3 - 1.5 mL		30B
HTRF Mouse IFN g kit - Std		05RA
HTRF Mouse IFN g kit - 500 pts d2 antibody		05RB
HTRF Mouse IFN g kit - 500 pts Eu Cryptate antibody		05RB

Results :

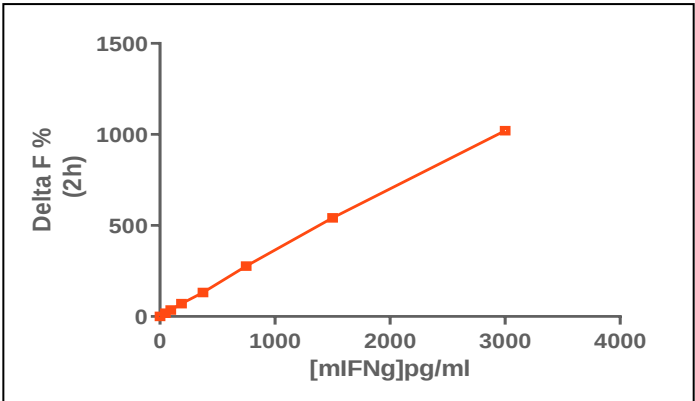
Incubation : 2h at RT

Plate : 384 wells small volume white plate 20 µL

Reader : Pherastar Plus

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

M IFN gamma	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	46,88	18%	13-24
Std 2	93,75	35%	26-42
Std 3	187,5	71%	62-75
Std 4	375	132%	114-168
Std 5	750	277%	220-358
Std 6	1500	542%	435-720
Std 7	3000	1021%	905-1279



Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	11 pg/ml	≤ 46.88 pg/ml

HTRF Mouse IFN g kit - 500 pts - Lot n° : 05G

☒ 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.