

HTRF® Research Reagents

HTRF Mouse IFN beta kit - 500 pts

Product Number : 62MIFNBPEG Lot Number : 03F

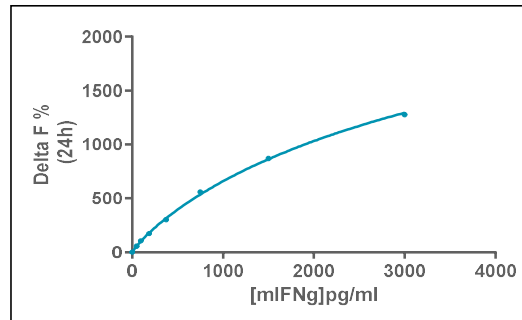
Expiration date : 2025-06-12 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21B
HTRF Detect. Buf. 3 - 1.5 mL	30B
HTRF Mouse IFN beta kit - Std	02A
HTRF Mouse IFN beta kit - 500 pts d2 antibody	03RA
HTRF Mouse IFN beta kit - 500 pts Eu Cryptate antibody	03RA

Results :

Incubation : Overnight at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Envision
 Fitting model : 4 Parameter Logistic (4PL 1/y2)* model

M IFN gamma	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	59,0	55%	50-76
Std 2	106	105%	85-128
Std 3	191	174%	156-235
Std 4	343	303%	267-401
Std 5	617	559%	459-689
Std 6	1111	871%	760-1140
Std 7	2000	1278%	1213-1820



Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	4,3 pg/ml	≤ 59 pg/ml

HTRF Mouse IFN beta kit - 500 pts - Lot n° : 03E

Accepted

Rejected

Quality control date : 2024-08-09	Supervisor's initials : MP
-----------------------------------	----------------------------

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.