

HTRF® Research Reagents

HTRF Human IFN beta kit - 10K pts

Product Number : 62HIFNBPEH Lot Number : 07A

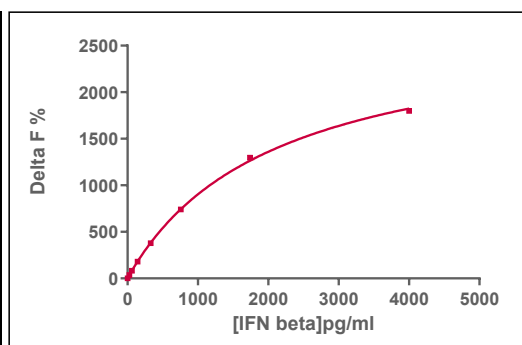
Expiration date : 2027-06-20 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 10 mL	22A
HTRF Activation reagent - 20 mL	04A
HTRF Detect. Buf. 3 - 50 mL	29RA
HTRF Human IFN beta kit - 10K pts d2 antibody	07A
HTRF Human IFN beta kit - 10K pts Eu Cryptate antibody	07A
HTRF Human IFN beta kit - Std	05RA
/	/

Results :

Incubation : 3h at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Nexus
 Fitting model : 4 Parameter Logistic (4PL 1/y2)* model

hIFN beta	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	27	37%	19--51%
Std 2	62	82%	54--115%
Std 3	143	181%	138--231%
Std 4	329	379%	273--493%
Std 5	756	741%	633--929%
Std 6	1739	1296%	1210--1579%
Std 7	4000	1798%	1773--2085%



	Found value	Specifications Rev 02
Detection Limit	2,9 pg/ ml	< 5.3pg/ml
	/	/

HTRF Human IFN beta kit - 10K pts - Lot n° : 07A

☒ Accepted

☐ Rejected

Quality control date : 2025-08-06	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.