

HTRF® Research Reagents

HTRF Human IFN beta kit - 10K pts

Product Number : 62HIFNBPEH Lot Number : 06A

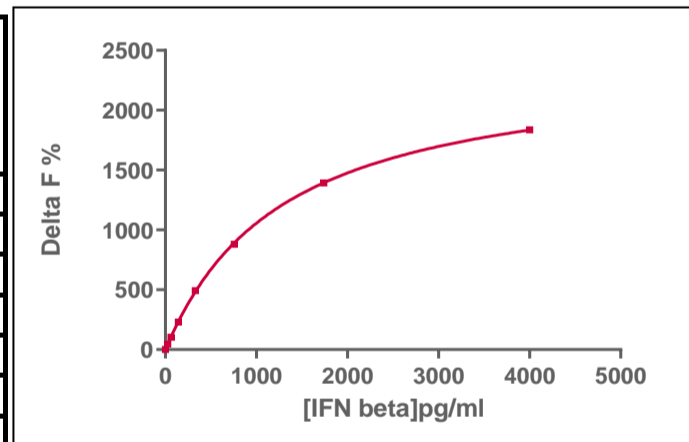
Expiration date : 2026-01-04 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 10 mL	21RA
HTRF Activation reagent - 20 mL	03RA
HTRF Detect. Buf. 3 - 50 mL	29RA
HTRF Human IFN beta kit - 10K pts d2 antibody	06A
HTRF Human IFN beta kit - 10K pts Eu Cryptate antibody	06A
HTRF Human IFN beta kit - Std	05RA

Results :

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

hIFN beta	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	27	46%	19--51%
Std 2	62	104%	54--115%
Std 3	143	230%	138--231%
Std 4	329	493%	273--493%
Std 5	756	881%	633--929%
Std 6	1739	1393%	1210--1579%
Std 7	4000	1837%	1773--2085%



	Found value	Specifications Rev 02
Detection Limit	<b>2,3 pg/ ml</b>	<b>&lt; 5.3pg/ml</b>
	/	/

HTRF Human IFN beta kit - 10K pts - Lot n° : 06A

Accepted

Rejected

Quality control date : 2024-08-23	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.