

HTRF® Research Reagents

### HTRF Human IFN beta kit - 500 pts

Product Number : 62HIFNBPEG Lot Number : 05J

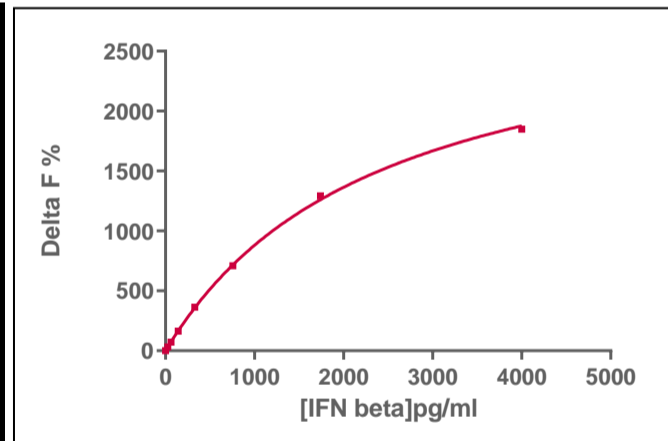
Expiration date : 2025-12-21 Document version : 1

Reagents :	Lot n°
HTRF Activation reagent - 1 mL	04A
HTRF Diluent 5 (5X) - 2 mL	21B
HTRF Detect. Buf. 3 - 1.5 mL	30B
HTRF Human IFN beta kit - Std	05RA
HTRF Human IFN beta kit - 500 pts d2 antibody	05RA
HTRF Human IFN beta kit - 500 pts Eu Cryptate antibody	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	40%	19--51%
Std 2	62	88%	54--115%
Std 3	143	188%	138--231%
Std 4	329	427%	273--493%
Std 5	756	821%	633--929%
Std 6	1739	1457%	1210--1579%
Std 7	4000	2065%	1773--2085%



**Assay characteristics :**

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

HTRF Human IFN beta kit - 500 pts - Lot n° : 05J

Accepted

Rejected

Quality control date : 2024-12-03	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.