

HTRF® Research Reagents

HTRF HP Human IL1 beta Kit - 96 pts

Product Number : 62HIL1B2PET Lot Number : 01B

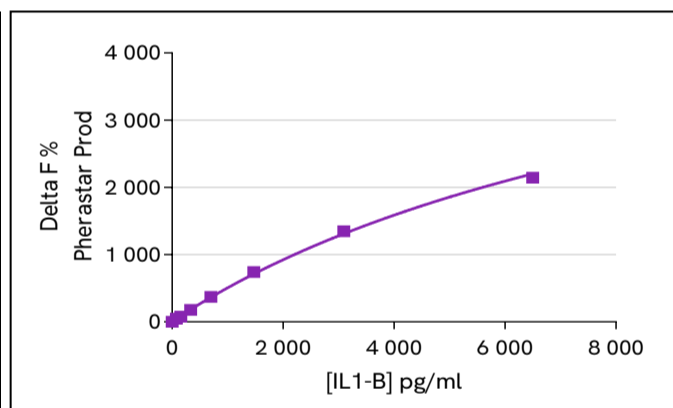
Expiration date : 2026-09-12 Document version : 1

Reagents :	Lot n°
PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND	2203
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 0.5 mL	30AR
HTRF HP Human IL1 beta Kit - Std	01A
HTRF HP Human IL1 beta Kit - 96 pts d2 antibody	01B
HTRF HP Human IL1 beta Kit - 96 pts Eu Cryptate antibody	01B

Results :

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

HP human IL1beta	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	75,8	48%	22-62
Std 2	159,1	78%	61-112
Std 3	334,2	177%	144-211
Std 4	701,9	370%	305-429
Std 5	1473,9	741%	605-891
Std 6	3095,2	1349%	1149-835
Std 7	6500	2144%	1722-2757



Assay characteristics :

	Found value	Specifications Rev 01
Detection limite	10 pg/ml	≤ 75,8pg/ml

HTRF HP Human IL1 beta Kit - 96 pts - Lot n° : 01B

Accepted

Rejected

Quality control date : 2024-07-03	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.