

HTRF® Research Reagents

HTRF HP Human IL1 beta Kit - 10K pts

Product Number : 62HIL1B2PEH Lot Number : 01A

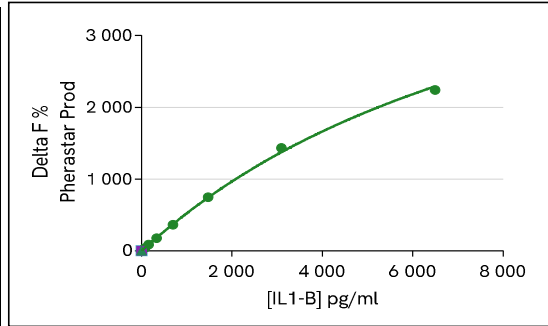
Expiration date : 2027-06-20 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 10 mL	21RA
HTRF Detect. Buf. 3 - 50 mL	29RA
HTRF HP Human IL1 beta Kit - 10K pts d2 antibody	01A
HTRF HP Human IL1 beta Kit - 10K pts Eu Cryptate antibody	01A
HTRF HP Human IL1 beta Kit - Std	01A
HTRF HP Human IL1 beta Kit - modop	2

Results :

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

HP human IL1beta	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	75.8	42%	22-62
Std 2	159.1	87%	61-112
Std 3	334.2	178%	144-211
Std 4	701.9	367%	305-429
Std 5	1473.9	748%	605-891
Std 6	3095.2	1434%	1149-835
Std 7	6500	2240%	1722-2757



Assay characteristics :

	Found value	Specifications Rev 01
Detection limite	pg/ml	≤ 159.1pg/ml

HTRF HP Human IL1 beta Kit - 10K pts - Lot n° : 01A

Accepted

Rejected

Quality control date : 2023-11-09	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.