

HTRF® Research Reagents

HTRF Human IL1 beta kit - 96 pts

Product Number : 62HIL1BPET Lot Number : 01ZB

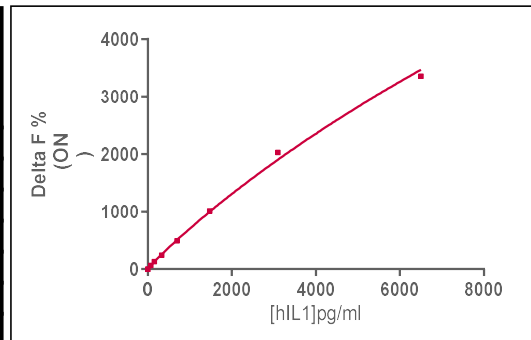
Expiration date : 2025-08-10 Document version : 1

Reagents :	Lot n°
PLAQUE BLANCHE 96 PUIITS SV COSMOBRAND	2413
HTRF Diluent 5 (5X) - 2 mL	21B
HTRF Detect. Buf. 3 - 0.5 mL	30AR
HTRF Human IL1 beta kit - Std	05RA
HTRF Human IL1 beta kit - 96 pts Eu Cryptate antibody	01RA
HTRF Human IL1 beta kit - 96 pts XL665 antibody	01RA

Results :

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

IL1 beta	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	75,8	66%	39-102
Std 2	159,1	130%	80-219
Std 3	334,2	244%	175-403
Std 4	701,9	495%	382-793
Std 5	1474	1010%	837-1508
Std 6	3095	2031%	1704-2800
Std 7	6500	3354%	2840-4975



Assay characteristics :

	Found value	Specifications Rev 02
Detection limite	4.5 pg/ml	≤ 75.8 pg/ml

HTRF Human IL1 beta kit - 96 pts - Lot n° : 01ZB

Accepted

Rejected

Quality control date :	2024-08-05	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 TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.