

HTRF® Research Reagents

HTRF Mouse IL1 beta kit - 500 pts

Product Number : 62MIL1BPEG Lot Number : 06F

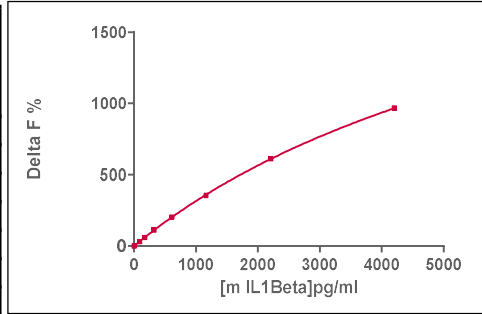
Expiration date : 2025-10-20 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21B
HTRF Detect. Buf. 3 - 1.5 mL	30B
HTRF Mouse IL1 beta kit - Std	05RA
HTRF Mouse IL1 beta kit - 500 pts d2 antibody	06RA
HTRF Mouse IL1 beta kit - 500 pts Eu Cryptate antibody	06RA

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

mIL1 beta	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	89	32%	31 - 47
Std 2	170	60%	52 - 91
Std 3	322	113%	103 - 162
Std 4	612	204%	186 - 297
Std 5	1163	356%	325 - 516
Std 6	2211	612%	549 - 821
Std 7	4200	968%	821 - 1067



Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	11 pg/ml	≤ 89 pg/ml

HTRF Mouse IL1 beta kit - 500 pts - Lot n° : 06F

Accepted

Rejected

Quality control date : 2027-07-02	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.