

HTRF® Research Reagents

HTRF Mouse IL1 beta kit - 10K pts

Product Number : 62MIL1BPEH Lot Number : 07A

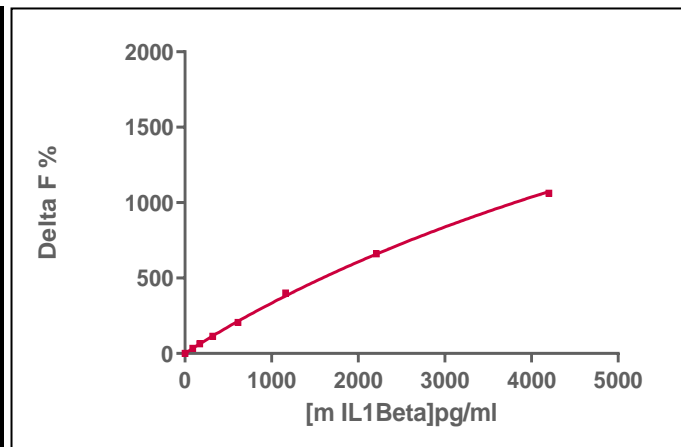
Expiration date : 2026-01-17 Document version : 1

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 10 mL	21RA
HTRF Detect. Buf. 3 - 50 mL	29RA
HTRF Mouse IL1 beta kit - 10K pts d2 antibody	07A
HTRF Mouse IL1 beta kit - 10K pts Eu Cryptate antibody	07A
HTRF Mouse IL1 beta kit - Std	05RA

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

m IL1 beta	Conc. (pg/ml)	DF %	DF Accepted Range
Std 1	89	34%	31 - 47
Std 2	170	64%	52 - 91
Std 3	322	114%	103 - 162
Std 4	612	205%	186 - 297
Std 5	1163	400%	325 - 516
Std 6	2211	661%	549 - 821
Std 7	4200	1063%	821 - 1067



Assay characteristics :

	Found value	Specifications Rev 01
Detection limite	10,6 pg/ml	≤ 89 pg/ml

HTRF Mouse IL1 beta kit - 10K pts - Lot n° : 07A

Accepted

Rejected

Quality control date : 2024-04-26	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.