

HTRF® Research Reagents

HTRF Mouse IL1 beta kit - 10K pts

Product Number : 62MIL1BPEH Lot Number : 06E

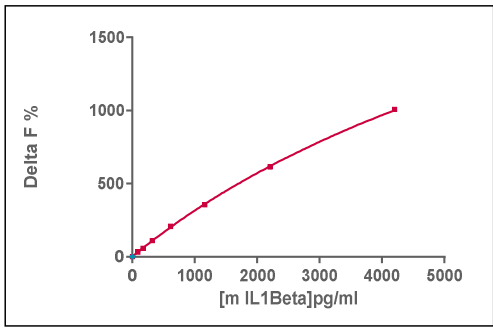
Expiration date : 2025/10/20 Document version : 1

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	06RA
HTRF Mouse IL1 beta kit - 10K pts Eu Cryptate antibody	06RA
HTRF Mouse IL1 beta kit - Std	05RA
HTRF Mouse IL1 beta kit - modop	7

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	32%	31 - 47
Std 2	170	58%	52 - 91
Std 3	322	111%	103 - 162
Std 4	612	208%	186 - 297
Std 5	1163	355%	325 - 516
Std 6	2211	614%	549 - 821
Std 7	4200	1008%	821 - 1067



Assay characteristics :

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

HTRF Mouse IL1 beta kit - 10K pts - Lot n° : 06E

Accepted

Rejected

Quality control date : 2023-09-08	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.