

HTRF® Research Reagents

HTRF (h) MMP1 Kit - 500 pts

Product Number : 62MMP1PEG Lot Number : 01B

Expiration date : 2026-10-16 Document version : 1

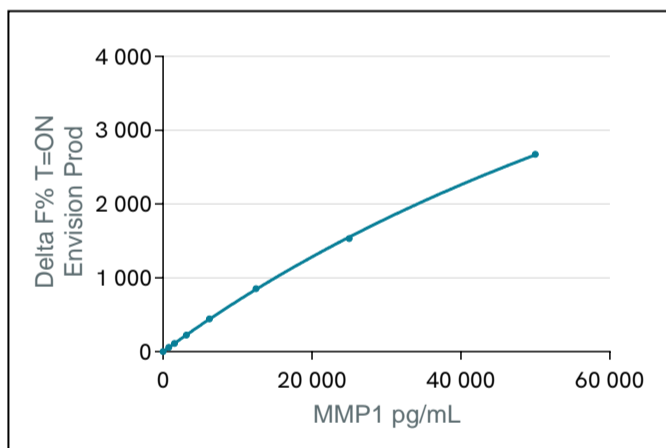
Reagents :		Lot n°
HTRF Diluent 5 (5X) - 2 mL		21B
HTRF Detect. Buf. 3 - 1.5 mL		30B
HTRF (h) MMP1 Kit - Std		01A
HTRF (h) MMP1 Kit - 500 pts d2 antibody		01A
HTRF (h) MMP1 Kit - 500 pts Eu Cryptate antibody		01A

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

Assay characteristics :

MMP1	Conc. (pg/ml)	DF %	DF Accepted Range
Std 0	0	0	0
Std 1	781	57	51-77
Std 2	1 563	113	102 - 153
Std 3	3 125	223	201 - 301
Std 4	6 250	443	382 - 573
Std 5	12 500	853	707 - 1061
Std 6	25000	1533	1241 - 1862
Std 7	50 000	2675	2088 - 3132



HTRF (h) MMP1 Kit - 500 pts - Lot n° : 01B

Accepted

Rejected

Quality control date : 2024-12-06	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.