

HTRF® Research Reagents

HTRF (h) MMP1 Kit - 500 pts

 Product Number : **62MMP1PEG** Lot Number : **01A**

 Expiration date : **16/10/2026** Document version : **1**

Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 1.5 mL	30AR
HTRF (h) MMP1 Kit - Std	01A
HTRF (h) MMP1 Kit - 500 pts d2 antibody	01A
2026	01A

Results :

Incubation :

Plate :

 Reader : **Pherastar Plus**

Fitting model :

Assay characteristics :

	Std en Pg/mL	Delta F (%)	Normes mini.	Normes maxii.
Std 0	0	0	0%	0%
Std 1	781	78	60%	96%
Std 2	1 563	157	123%	192%
Std 3	3 125	317	277%	357%
Std 4	6 250	602	550%	655%
Std 5	12 500	1110	1071%	1150%
Std 6	25000	1982	1923%	2041%
Std 7	50 000	3274	2940%	3607%

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

 Rejected

Quality control date : 2024-02-12	Supervisor's initials : AC
--	-----------------------------------

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.