

## **MMP-1 Colorimetric Drug Discovery Kit**

Catalog #: [BML-AK404](#)

Designed to screen MMP-1 inhibitors using a thiopeptide as a chromogenic substrate.

## **USE FOR RESEARCH PURPOSES ONLY**

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

## **LIMITED WARRANTY; DISCLAIMER OF WARRANTIES**

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NON-INFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

## **TRADEMARKS AND PATENTS**

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending. Enzo is a trademark of Enzo Life Sciences, Inc.

**FOR RESEARCH USE ONLY.  
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**



Please read entire booklet before proceeding with the assay.



Carefully note the handling and storage conditions of each kit component.



Please contact Enzo Life Sciences Technical Support if necessary.

## TABLE OF CONTENTS

Background .....	2
References .....	2
Materials Supplied, Storage .....	3
Other Materials Needed .....	4
Experimental Methods.....	5
Data Analysis .....	7
Contact Information .....	10

## BACKGROUND

Matrix metalloproteinase-1 (MMP-1, interstitial collagenase, fibroblast collagenase) is a member of the MMP family of extracellular proteases. These enzymes play a role in many normal and disease states by virtue of their broad substrate specificities<sup>1</sup>. Targets of MMP-1 include collagen, gelatin, entactin, pro-TNF- $\alpha$ , and the chemokine SDF-11-4. MMP-1 is secreted as a 52-56kDa glycosylated proenzyme (as measured by SDS-PAGE), and activated by cleavage to forms of 22-46kDa<sup>5</sup>. MMP-1 is an important target for inhibitor screening due to its involvement in diseases such as cancer<sup>3</sup>.

The *MMP-1 Colorimetric Drug Discovery Kit* is a complete assay system designed to screen MMP-1 inhibitors using a thiopeptide as a chromogenic substrate (Ac-PLG-[2-mercapto-4-methyl-pentanoyl]-LG-OC<sub>2</sub>H<sub>5</sub>)<sup>6,7</sup>. The MMP cleavage site peptide bond is replaced by a thioester bond in the thiopeptide. Hydrolysis of this bond by an MMP produces a sulfhydryl group, which reacts with DTNB [5,5'-dithiobis(2-nitrobenzoic acid), Ellman's reagent] to form 2-nitro-5-thiobenzoic acid, which can be detected by its absorbance at 412 nm ( $\epsilon=13,600 \text{ M}^{-1}\text{cm}^{-1}$  at pH 6.0 and above<sup>8</sup>). The assays are performed in a convenient 96-well microplate format. The kit is useful to screen inhibitors of MMP-1, a potential therapeutic target. An inhibitor, NNGH9, is also included as a prototypic control inhibitor. Thiol inhibitors should not be used with this kit, as they may interfere with the colorimetric assay.

## REFERENCES:

1. L.J. McCawley and L.M. Matrisian Curr. Opin. Cell Biol. 2001 13 534
2. H. Nagase and J.F. Woessner, Jr. J. Biol. Chem. 1999 274 2149
3. W.G. Stetler-Stevenson and A.E. Yu Semin. Cancer Biol. 2001 11 143
4. G.A. McQuibban et al. J. Biol. Chem. 2001 276 43503
5. K. Suzuki et al. Biochemistry 1990 29 10261
6. H. Weingarten and J. Feder Anal. Biochem. 1985 147 437
7. H. Weingarten et al. Biochemistry 1985 24 6730
8. L. Yu and E.A. Dennis Methods Enzymol. 1991 197 65
9. L.J. MacPherson et al. J. Med. Chem. 1997 40 2525



Avoid  
freeze /  
thaw cycles



Reagents  
require  
separate  
storage  
conditions.

## MATERIALS SUPPLIED

- 1. MMP-1 ENZYME (HUMAN, RECOMBINANT)**  
**Catalog No. BML-SE180-9090:**  
**FORM:** E. coli recombinant human MMP-1 catalytic domain (calculated MW 19.9kDa), 30.6 U/ $\mu$ l. One U=100 pmol/min@ 37°C, 100 $\mu$ M thiopeptide P125.  
**STORAGE:** -70°C; Avoid freeze/thaw cycles  
**QUANTITY:** 2000 U
- 2. MMP SUBSTRATE (chromogenic; MW=655.9)**  
**Catalog No. BML-P125-9090:**  
**FORM:** 25mM (16.4mg/ml) in DMSO  
**STORAGE:** -20°C  
**QUANTITY:** 50 $\mu$ l
- 3. INHIBITOR (NNGH; MW=316.4)**  
**Catalog No. BML-PI115-9090:**  
**FORM:** 1.3mM in DMSO  
**STORAGE:** -20°C  
**QUANTITY:** 50 $\mu$ l
- 4. ASSAY BUFFER, Catalog No. BML-KI173-0020:**  
50mM HEPES, 10mM CaCl<sub>2</sub>, 0.05% Brij-35, 1mM DTNB, pH7.5  
**STORAGE:** -20°C  
**QUANTITY:** 20 ml
- 5. 96-WELL MICROPLATE, Catalog No. 80-2404:**  
**STORAGE:** Room temperature.

## **OTHER MATERIALS NEEDED**

1. Microplate reader capable of measuring A412 to  $\geq 3$ -decimal accuracy.
2. Pipetman or multi-channel pipetman capable of pipetting 10-100 $\mu$ l accurately (note: reagents can be diluted to increase the minimal pipetting volume to  $>10\mu$ l).
3. Ice bucket to keep reagents cold until use.
4. Water bath or incubator for component temperature equilibration.

## EXPERIMENTAL METHODS

**Note on storage:** Store all components except the microplate (room temperature) at -70°C for the highest stability. The MMP-1 enzyme should be handled carefully in order to retain maximal enzymatic activity. It is stable, in diluted or concentrated form, for several hours on ice. As supplied, MMP-1 enzyme is stable for at least 4 freeze/thaw cycles. To minimize the number of freeze/thaw cycles, aliquot the MMP-1 into separate tubes and store at -70°C. When setting up the assay, do not maintain diluted components at reaction temperature (e.g., 37°C) for an extended period of time prior to running the assay.

### To start assay:

1. Briefly warm kit components BML-P125-9090 and BML-PI115-9090 to RT to thaw DMSO.
2. Dilute inhibitor (NNGH, BML-PI115-9090) 1/200 in assay buffer BML-KI173 as follows. Add 1µl inhibitor into 200µl assay buffer, in a separate tube. Warm to reaction temperature (e.g., 37°C).
3. Dilute substrate BML-P125-9090 1/25 in assay buffer to required total volume (10µl are needed per well). For example, for 15 wells dilute 6.4µl BML-P125-9090 into 153.6µl assay buffer, in a separate tube. Warm to reaction temperature (e.g., 37°C).
4. Dilute MMP-1 enzyme 1/40 in assay buffer to required total volume (20µl are needed per well). Warm to reaction temperature (e.g., 37°C) shortly before assay.
5. Pipet assay buffer into each desired well of the 1/2 volume microplate as follows:  
**Blank** (no MMP-1)=90µl Assay Buffer  
**Control** (no inhibitor)=70µl Assay Buffer  
**Inhibitor** NNGH=50µl Assay Buffer  
**Test inhibitor**=varies (see Table 1, below)
6. Allow microplate to equilibrate to assay temperature (e.g., 37°C).
7. Add 20µl MMP-1 (diluted in step 4) to the control, inhibitor NNGH, and test inhibitor wells. Final amount of MMP-1 will be 15.3 U per well (153 mU/µl). Remember to not add MMP-1 to the blanks!
8. Add 20µl NNGH inhibitor (diluted in step 2) to the inhibitor NNGH wells only! Final inhibitor concentration=1.3µM. **Note:** 1µM NNGH will not completely inhibit MMP-1 under these conditions (see **Figure 2**).
9. Add desired volume of test inhibitor to appropriate wells. See **Table 1**, following.
10. Incubate plate for 30-60 minutes at reaction temperature (e.g., 37°C) to allow inhibitor/enzyme interaction.

11. Start reaction by the addition of 10µl BML-P125-9090 substrate (diluted and equilibrated to reaction temperature in step 3). Final substrate concentration=100µM.
12. Continuously read plates at A412nm in a microplate reader. Record data at 1 min. time intervals for 10 to 20 min.
13. Perform data analysis (see below).

**NOTE:** Retain microplate for future use of unused wells.

**Table 1. Example of Samples.**

Sample	Assay buffer	MMP-1 (765mU/µl)	Inhibitor (6.5µM)	Substrate (1mM)	Total Volume
Blank	90µl	0	0	10µl	100µl
Control	70µl	20µl	0	10µl	100µl
Inhibitor NNGH	50µl	20µl	20µl	10µl	100µl
Test inhibitor*	Xµl	20µl	Yµl	10µl	100µl

\*Test inhibitor is the experimental inhibitor. Dissolve/dilute inhibitor into assay buffer and add to appropriate wells at desired volume “Y”. Adjust volume “X” to bring the total volume to 100µl.

Example of plate:	<u>well#</u>	<u>Sample</u>
	A1	Blank
	B1	Blank
	C1	Control
	D1	Control
	E1	Inhibitor NNGH
	F1	Inhibitor NNGH
	G1	Test inhibitor
	H1	Test inhibitor



## DATA ANALYSIS

### Plotting

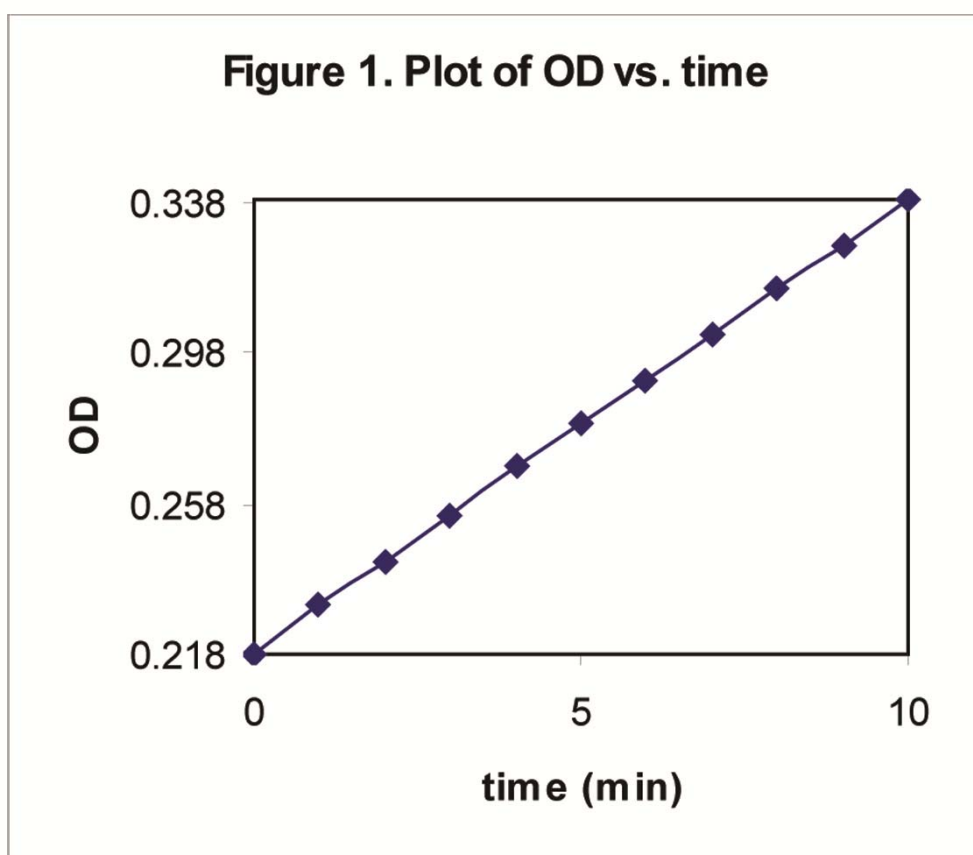
1. Plot data as OD versus time for each sample (see **Fig. 1**).
2. Determine the range of time points during which the reaction is linear. Typically, points from 1 to 10 min are sufficient.
3. Obtain the reaction velocity (V) in OD/min: determine the slope of a line fit to the linear portion of the data plot using an appropriate routine.
4. Average the slopes of duplicate samples.

### Data Reduction

5. If the blank has a significant slope, subtract this number from all samples.
6. To determine inhibitor % remaining activity:

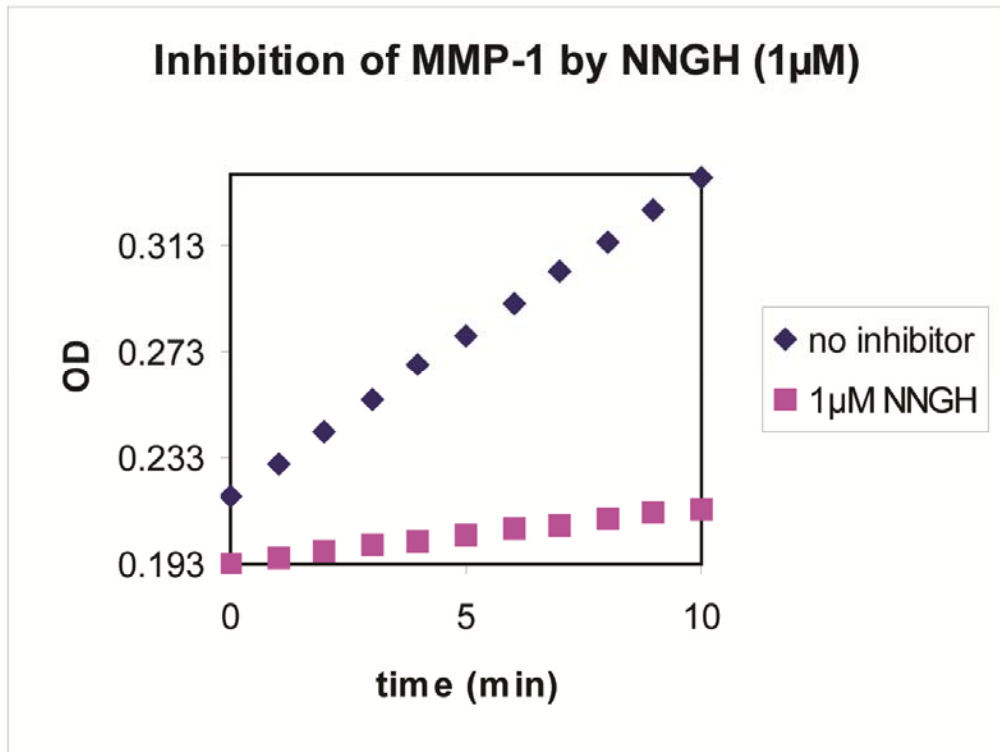
$$\text{Inhibitor \% activity remaining} = (\mathbf{V} \text{ inhibitor} / \mathbf{V} \text{ control}) \times 100$$

See **Figure 2** for example.



Slope=V=1.20E-02 OD/min

**Figure 2**

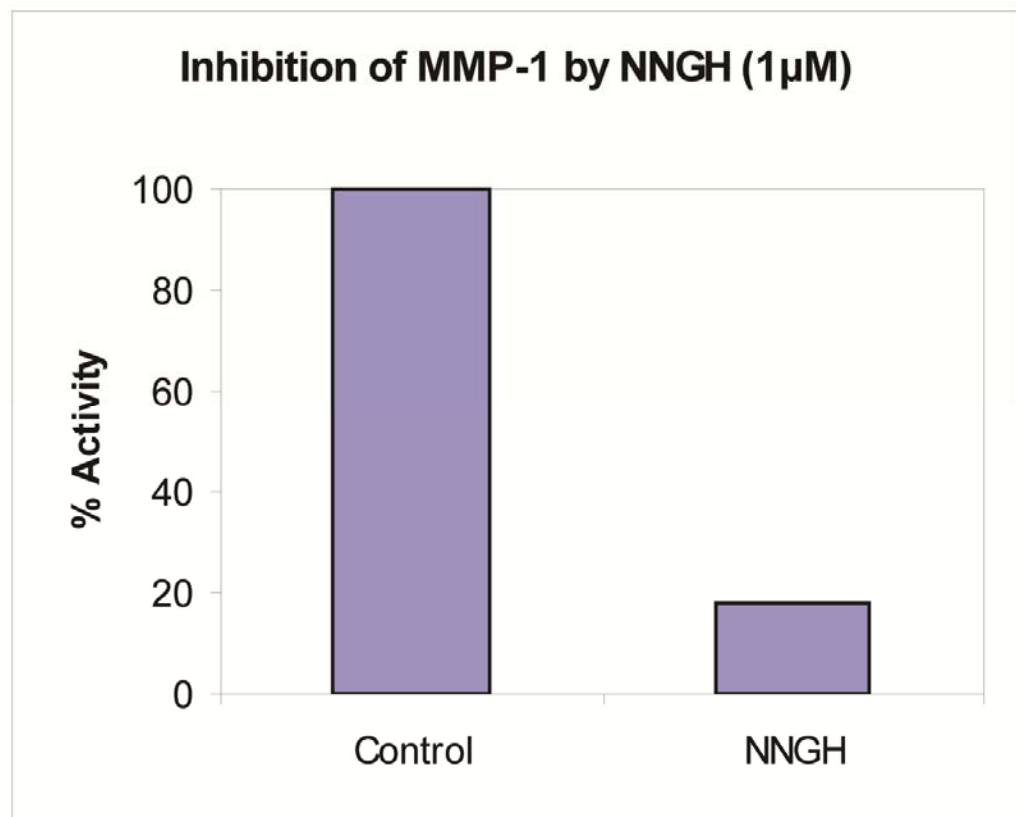


control slope = 1.20E-02 OD/min

inhibitor slope = 2.13E-03 OD/min

inhibitor % activity remaining =  $(2.13E-03/1.20E-02) \times 100 = 17.7\%$

**Figure 3**



7. To find the activity of the samples expressed as mol substrate/min, employ the following equation:

$$X \text{ mol substrate/min} = (V \times \text{vol.}) / (\epsilon \times \ell)$$

Where **V** is reaction velocity in OD/min, vol. is the reaction volume in liters,  $\epsilon$  is the extinction coefficient of the reaction product (2-nitro-5-thiobenzoic acid) ( $13,600 \text{ M}^{-1}\text{cm}^{-1}$ ), and  $\ell$  is the path length of light through the sample in cm (for  $100\mu\text{l}$  in the supplied microplate,  $\ell$  is 0.5 cm).

**Note:** The above equation determines enzyme activity in terms of moles of thiopeptolide substrate P125 converted per minute. Under these conditions, the secondary substrate DTNB is saturating, and the velocity of DTNB conversion to 2-nitro-5-thiobenzoic acid is not rate-limiting.

See **Figure 3** for activity and kinetic calculations.

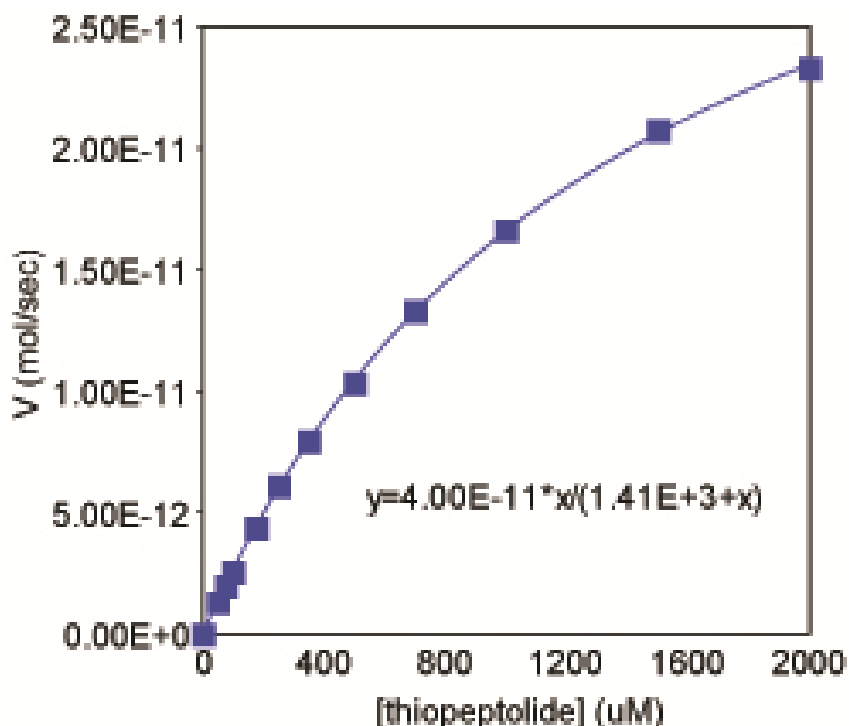
**Figure 3. Example calculation for activity:**

Activity of a control sample =

$$(1.04\text{E-}02\text{OD/min} \times 1\text{E-}04\text{L}) / (13,600\text{M}^{-1}\text{cm}^{-1} \times 0.5\text{cm}) = 1.53\text{E-}10 \text{ mol/min at } 37^\circ\text{C, } 100\mu\text{M thiopeptolide P125}$$

**Example graph for  $K_m$  and  $V_{max}$  determination:**

**Figure 4**



$$K_m = 1410\mu\text{M}$$

$$V_{max} = 40.0 \text{ pmol/sec}$$

**GLOBAL HEADQUARTERS**

Enzo Life Sciences Inc.  
10 Executive Boulevard  
Farmingdale, NY 11735  
Toll-Free: 1.800.942.0430  
Phone: 631.694.7070  
Fax: 631.694.7501  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

**EUROPE/ASIA**

Enzo Life Sciences (ELS) AG  
Industriestrasse 17  
CH-4415 Lausen  
Switzerland  
Phone: +41/0 61 926 89 89  
Fax: +41/0 61 926 89 79  
[info-ch@enzolifesciences.com](mailto:info-ch@enzolifesciences.com)

For local distributors and detailed product information visit us online:  
[www.enzolifesciences.com](http://www.enzolifesciences.com)