



Product Manual

MMP Inhibitor Profiling Kit, Colorimetric

Designed to examine the specificity of inhibitors against a panel of ten matrix metalloproteinase enzymes.

Instruction Manual

BML-AK015

For research use only



Product Manual

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.**

BACKGROUND

The matrix metalloproteinases, or MMPs, are extracellular proteases that function at a neutral pH to cleave a wide variety of substrates. These include basement membrane and extracellular matrix components, growth and death factors, cytokines, and cell and matrix adhesion molecules^{1,2,3}. The broad range of substrate specificities and expression patterns of MMPs results in their involvement in many different processes, both normal and pathological. Aberrant expression has been noted in cancer, angiogenesis, arthritis, inflammation, periodontal disease, emphysema, multiple sclerosis, pre-eclampsia, and chronic wounds, among others^{1,2,4}. The general structure of an MMP protein consists of a pro domain to direct secretion from the cell, a catalytic domain, and a C-terminal hemopexin domain. The catalytic site involves a coordinately-bound zinc ion. The inactive, or zymogen, form of the enzyme is activated by disruption of one of the coordinate bonds, usually via proteolytic removal of the pro domain⁵.

The *MMP Inhibitor Profiling Kit* is a complete assay system designed to examine the specificity of inhibitors against a panel of ten matrix metalloproteinase enzymes. A thiopeptide is used as a chromogenic substrate (Ac-PLG-[2-mercapto-4-methylpentanoyl]-LG-OC₂H₅)^{6,7}. 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 above⁸). The assays are performed in a convenient 96-well microplate format. 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.

Please contact Enzo Life Sciences for kit components in bulk.

REFERENCES:

1. G. Bergers and L.M. Coussens *Curr. Opin. Genet. Dev.* 2000 **10** 120
2. S.D. Shapiro *Curr. Opin. Cell Biol.* 1998 **10** 602
3. V. Noe *et al. J. Cell Sci.* 2001 **114** 111
4. A.R. Nelson *et al. J. Clin. Oncol.* 2000 **18** 1135
5. J.F. Woessner and H. Nagase *Metalloproteinases and TIMPs.* 2000 Oxford University Press
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

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 SERVICES FOR ASSISTANCE IF NECESSARY.

MATERIALS SUPPLIED

BML-SE180-9090 MMP-1 ENZYME (HUMAN, RECOMBINANT)

FORM: Recombinant human MMP-1 catalytic domain (19.9kDa); 30.6U/ μ L. One U=100 pmol/min@ 37°C, 100 μ M thiopeptide P125.

STORAGE: -80°C; Avoid freeze/thaw cycles

QUANTITY: 2000U

BML-SE237-9090 MMP-2 ENZYME (HUMAN, RECOMBINANT)

FORM: Recombinant human MMP-2 catalytic domain (40kDa); 3.28U/ μ L. One U=100 pmol/min@ 37°C, 100 μ M thiopeptide.

STORAGE: -80°C; Avoid freeze/thaw cycles

QUANTITY: 150U

BML-SE109-9090 MMP-3 ENZYME (HUMAN, RECOMBINANT)

FORM: Recombinant human MMP-3 catalytic domain (19.5kDa); 10U/ μ L. One U=100 pmol/min@ 37°C, 100 μ M thiopeptide.

STORAGE: -80°C; Avoid freeze/thaw cycles

QUANTITY: 300U

BML-SE181-9090 MMP-7 ENZYME (HUMAN, RECOMBINANT)

FORM: Recombinant human MMP-7 catalytic domain (20.4kDa); 4.5U/ μ L. One U=100 pmol/min@ 37°C, 100 μ M thiopeptide.

STORAGE: -80°C; Avoid freeze/thaw cycles

QUANTITY: 150U

BML-SE255-9090 MMP-8 ENZYME (HUMAN, RECOMBINANT)

FORM: Recombinant human MMP-8 catalytic domain (20.3kDa); 9.2U/ μ L. One U=100 pmol/min@ 37°C, 100 μ M thiopeptide.

STORAGE: -80°C; Avoid freeze/thaw cycles

QUANTITY: 200U

BML-SE360-9090 MMP-9 ENZYME (HUMAN, RECOMBINANT)

FORM: Recombinant human MMP-9 catalytic domain (39kDa); 2.68U/ μ L. One U=100 pmol/min@ 37°C, 100 μ M thiopeptide.

STORAGE: -80°C; Avoid freeze/thaw cycles

QUANTITY: 130U

BML-SE329-9090 MMP-10 ENZYME (HUMAN, RECOMBINANT)

FORM: Recombinant human MMP-10 catalytic domain (19.4kDa); 5U/ μ L. One U=100 pmol/min@ 37°C, 100 μ M thiopeptide.

STORAGE: -80°C; Avoid freeze/thaw cycles

QUANTITY: 150U

BML-SE138-9090 MMP-12 ENZYME (HUMAN, RECOMBINANT)

FORM: Recombinant human MMP-12 catalytic domain (20.3kDa); 10U/ μ L. One U=100 pmol/min@ 37°C, 100 μ M thiopeptide.

STORAGE: -80°C; Avoid freeze/thaw cycles

QUANTITY: 140U

BML-SE246-9090 MMP-13 ENZYME (HUMAN, RECOMBINANT)

FORM: Recombinant human MMP-13 catalytic domain (20.4kDa); 3.45U/ μ L. One U=100 pmol/min@ 37°C, 100 μ M thiopeptide.

STORAGE: -80°C; Avoid freeze/thaw cycles

QUANTITY: 180U

BML-SE259-9090 MMP-14 ENZYME (HUMAN, RECOMBINANT)

FORM: Recombinant human MMP-14 catalytic domain (22.5kDa); 12U/ μ L. One U=100 pmol/min@ 37°C, 100 μ M thiopeptide.

STORAGE: -80°C; Avoid freeze/thaw cycles

QUANTITY: 300U

BML-P125-9090 MMP SUBSTRATE (chromogenic; MW=655.9)

FORM: 25mM (16.4mg/ml) in DMSO

STORAGE: -20°C

QUANTITY: 50 μ L

BML-PI115-9090 INHIBITOR (NNGH; MW=316.4)

FORM: 1.3mM in DMSO

STORAGE: -20°C

QUANTITY: 50 μ L

BML-KI173-0020 ASSAY BUFFER

FORM: 50mM HEPES, 10mM CaCl₂, 0.05% Brij-35, 1mM DTNB, pH7.5

STORAGE: -20°C

QUANTITY: 20 ml

80-2404 ½ Volume clear 96-Well Microplate

STORAGE: Ambient

OTHER MATERIALS REQUIRED

1. Microplate reader capable of measuring A_{412} to ≥ 3 -decimal accuracy.
2. Pipetman or multi-channel pipetman capable of pipetting 10-100 μ 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 at -80°C for the highest stability. Components with storage temperatures other than -80°C can be stored at the temperature listed OR at -80°C . The MMP enzymes should be handled carefully in order to retain maximal enzymatic activity. They are stable, in diluted or concentrated form, for several hours on ice. As supplied, the MMP enzymes are stable for at least 5 freeze/thaw cycles. If necessary, to minimize the number of freeze/thaw cycles, aliquot the MMPs into separate tubes and store at -80°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 P125-9090 and P1115-9090 to RT to thaw DMSO.
2. Dilute control inhibitor (NNGH, P1115-9090) 1/200 in assay buffer K1173 as follows. Add 1 μ L inhibitor into 200 μ L assay buffer, in a separate tube. Warm to reaction temperature (e.g., 37°C).
3. Dilute substrate P125-9090 1/25 in assay buffer to required total volume (10 μ L are needed per well). For example, for 15 wells dilute 6.4 μ L P125-9090 into 153.6 μ L assay buffer, in a separate tube. Warm to reaction temperature (e.g., 37°C).

4. Shortly before assay, dilute MMP enzymes in assay buffer to required total volume ($20\mu\text{L}$ are needed per well), using the following ratios for dilution:

MMP-1: 1/80	MMP-9: 1/60
MMP-2: 1/70	MMP-10: 1/100
MMP-3: 1/70	MMP-12: 1/285
MMP-7: 1/70	MMP-13: 1/50
MMP-8: 1/100	MMP-14: 1/100

Note: MMP-3 is unique in that its pH optimum is 6.0. Activity of this enzyme in pH 7.5 buffer is somewhat reduced; thus the dilution for MMP-3 has been adjusted accordingly.

Warm to reaction temperature (e.g., 37°C).

5. Pipet assay buffer into each desired well of the $1/2$ volume microplate as follows:

Blank (no MMP)= $90\mu\text{L}$ Assay Buffer

Control (no inhibitor)= $70\mu\text{L}$ Assay Buffer

Inhibitor NNGH= $50\mu\text{L}$ Assay Buffer

Test inhibitor=varies (see Table 1, below)

6. Allow microplate to equilibrate to assay temperature (e.g., 37°C).
7. Add $20\mu\text{L}$ MMP (diluted in step 4) to the control, inhibitor NNGH, and test inhibitor wells.
8. Add $20\mu\text{L}$ NNGH inhibitor (diluted in step 2) to the inhibitor NNGH wells only! Final inhibitor concentration= $1.3\mu\text{M}$. See Figure 2 for example of inhibition of MMPs by NNGH.
9. Add desired volume of test inhibitor to appropriate wells. See Table 1, below.
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 $10\mu\text{L}$ P125-9090 substrate (diluted and equilibrated to reaction temperature in step 3). Final substrate concentration= $100\mu\text{M}$.
12. Continuously read plates at $A_{412\text{nm}}$ 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	Inhibitor (6.5 μ M)	Substrate (1mM)	Total Volume
Blank	90 μ L	0	0	10 μ L	100 μ L
Control	70 μ L	20 μ L	0	10 μ L	100 μ L
Inhibitor NNGH	50 μ L	20 μ L	20 μ L	10 μ L	100 μ L
Test inhibitor*	X μ L	20 μ L	Y μ L	10 μ L	100 μ L

*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 100 μ L.

Example of plate:	<u>well#</u>	<u>sample</u>
	A1	Blank
	B1	Blank
	C1	Control/MMP-1
	D1	Control/MMP-1
	E1	Inhibitor NNGH/MMP-1
	F1	Inhibitor NNGH/MMP-1
	G1	Test inhibitor/MMP-1
	H1	Test inhibitor/MMP-1
	A2	Control/MMP-2
	B2	Control/MMP-2
	C2	Test inhibitor/MMP-2
	D2	Test inhibitor/MMP-2

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.

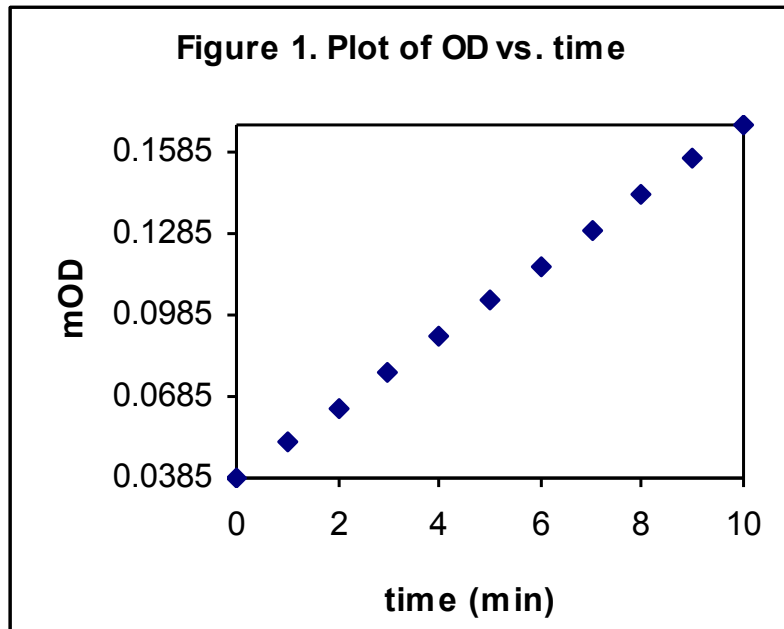
Data Reduction

5. *If the blank has a significant slope, subtract this number from all samples.*

6. To determine inhibitor % remaining activity:

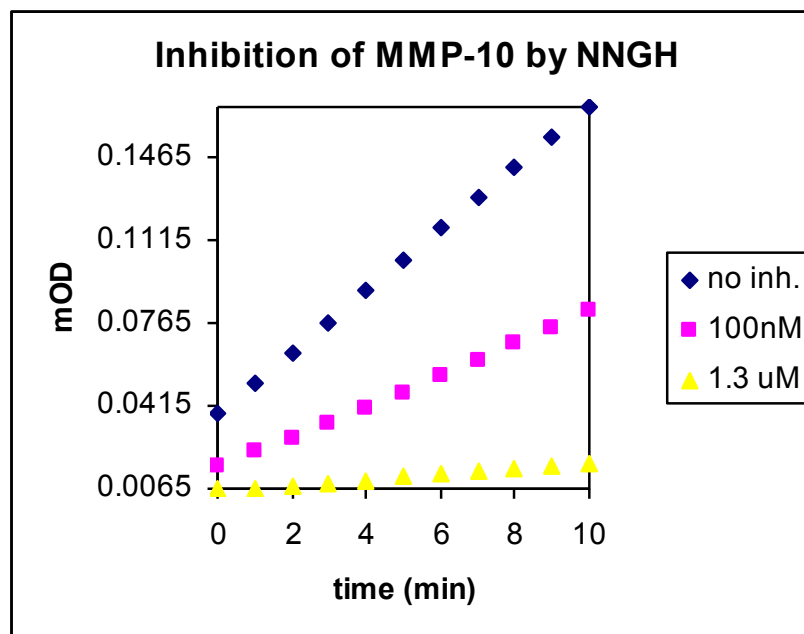
$$\text{Inhibitor \% activity remaining} = (\text{V inhibitor} / \text{V control}) \times 100$$

See **Figure 2** for example.



Slope=V=1.30E-02 OD/min

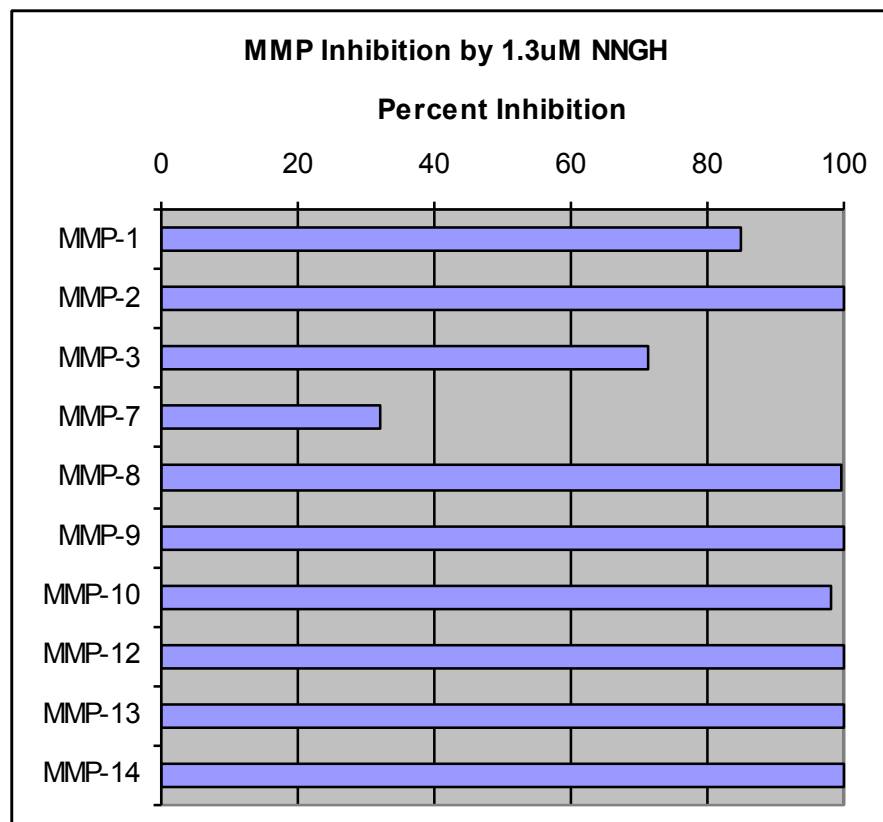
Figure 2. Example of inhibitor data:



control slope = 1.30E-02 OD/min

inhibitor (100nM) slope = 6.64E-03 OD/min

inhibitor % activity remaining = $(6.63\text{E-}03 / 1.30\text{E-}02) \times 100 = 51.0\%$



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, ϵ is the extinction coefficient of the reaction product (2-nitro-5-thiobenzoic acid) ($13,600 \text{ M}^{-1} \text{ cm}^{-1}$), and ℓ is the path length of light through the sample in cm (for 100ml in the supplied microplate, ℓ is 0.5 cm).

Note: The above equation determines enzyme activity in terms of moles of thiopeptide 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.

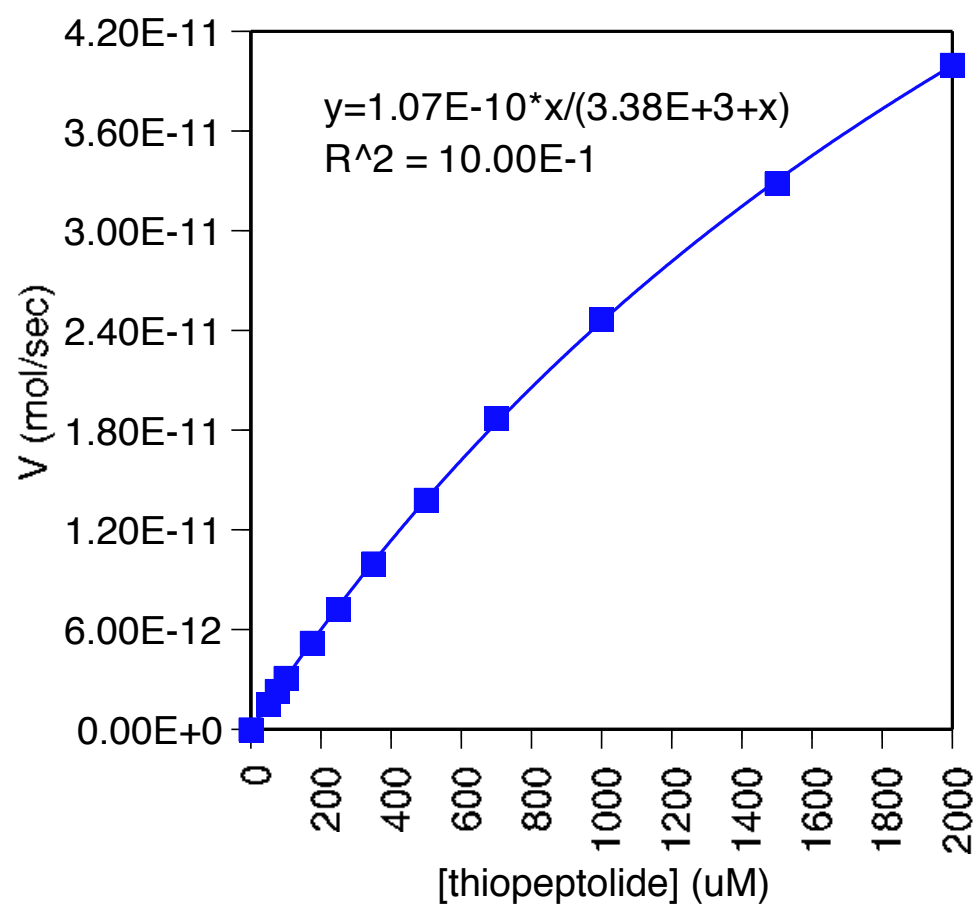
See **Figure 3** (following page) for activity and kinetic calculations.

Figure 3. Example calculation for activity:

Activity of a control sample =

$$(1.25E-02 \text{ OD/min} \times 1E-04 \text{ L}) / (13,600 \text{ M}^{-1} \text{ cm}^{-1} \times 0.5 \text{ cm}) = 1.84E-10 \text{ mol/L/min at } 37^\circ\text{C, } 100\mu\text{M thiopeptolide P125}$$

Example graph for K_m and V_{max} determination for MMP-10 with substrate P-125:



$K_m = 3380 \mu\text{M}$

$V_{max} = 107 \text{ pmol/sec}$



Product Manual

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

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

For local distributors and detailed product information visit us online:
www.enzolifesciences.com