



Enabling Discovery in Life Science®

MMP-3 Fluorometric Drug Discovery Kit

Designed to screen MMP-3 inhibitors using a quenched fluorogenic peptide.

**Instruction Manual
BML-AK401**

For research use only

BACKGROUND

Matrix metalloproteinase-3 (MMP-3, stromelysin-1, transin-1) 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 specificities¹. Targets of MMP-3 include collagens, fibronectin, and laminin, plasminogen, HB-EGF, E-cadherin, and other MMPs^{1,2,3}. MMP-3 is secreted as a 55-59 kDa glycosylated proenzyme (measured by SDS-PAGE), and activated by cleavage to forms of 21-48 kDa. It is unique from other MMPs in that its pH optimum is 5.9, rather than around 7.0⁴.

The *MMP-3 Fluorimetric Drug Discovery Kit* is a complete assay system designed to screen MMP-3 inhibitors using a quenched fluorogenic peptide: OmniMMP™ fluorogenic substrate Mca-Pro-Leu-Gly-Leu-Dpa-Ala-Arg-NH₂ [Mca=(7-methoxycoumarin-4-yl)-acetyl; Dpa=N-3-(2,4-dinitrophenyl)-L- α - β -diaminopropionyl]⁵. Mca fluorescence is quenched by the Dpa group until cleavage by MMPs at the Gly-Leu bond separates the two moieties^{5,6}. The assays are performed in a convenient 96-well microplate format. The kit is useful to screen inhibitors of MMP-3, a potential therapeutic target. The compound NNGH⁷ is also included as a prototypic control inhibitor.

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. Noë *et al. J. Cell Sci.* 2000 **114** 111
4. L.L. Johnson *et al. J. Biol. Chem.* 2000 **275** 11026
5. C.G. Knight *et al. FEBS Lett.* 1992 **296** 263
6. M.M. Smith *et al. J. Biol. Chem.* 1995 **270** 6440
7. 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.

COMPONENTS OF BML-AK401 KIT

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

FORM: *E. coli* recombinant human MMP-3 catalytic domain (calculated MW 19.5 kDa), 10 U/ μ l

UNIT DEFINITION: One unit is defined as the amount of enzyme that will hydrolyze 100 μ M thiopeptolide Ac-PLG-[2-mercapto-4-methyl-pentanoyl]-LG-OC₂H₅ (Cat. # BML-P125) at 100 pmol/min@ 37°C.

STORAGE: -70°C; Avoid freeze/thaw cycles

QUANTITY: 300 U

PRESENTATION: 30 μ l in screw-cap microfuge vial.

BML-P126-9090 SUBSTRATE (OmniMMP™ fluorogenic substrate peptide; MW=1093.2)

FORM: 400 μ M (437 μ g/ml) in DMSO (dimethylsulfoxide)

STORAGE: -70°C

PRESENTATION: 200 μ l in amber screw-cap microfuge vial.

BML-P127-9090 CALIBRATION STANDARD (OmniMMP™ fluorogenic control peptide, Mca-Pro-Leu-OH; MW=444.5)

FORM: 40 μ M (17.8 μ g/ml) in DMSO

STORAGE: -70°C

PRESENTATION: 50 μ l in amber screw-cap microfuge vial.

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

FORM: 1.3mM in DMSO

STORAGE: -20 or -70°C

PRESENTATION: 50 μ l in screw-cap microfuge vial.

BML-KI127-0020 ASSAY BUFFER

50mM MES, 10mM CaCl₂, 0.05% Brij-35, pH 6.0

FORM: Liquid in screw-cap plastic bottle

STORAGE: Room temperature

QUANTITY: 20 ml

80-2406 1/2 VOLUME 96-WELL WHITE NBS MICROPLATE

STORAGE: Room temperature.

(Sold separately as BML-KI571)

OTHER MATERIALS REQUIRED

Fluorescent microplate reader capable of excitation at 328nm and emission at 420nm. The following Ex/Em have also been used: 320,340/393,400,405.

Pipetmen or multi-channel pipetmen capable of pipetting 1-100 μ l accurately

Ice bucket to keep reagents cold until use.

Water bath or incubator for component temperature equilibration

EXPERIMENTAL METHODS

PLEASE NOTE: *The recommended dilution of substrate BML-P126-9090 in step 3 has been changed so that the volume added to assay wells (in steps 11 and 12) is 10 μ l instead of 1 μ l.*

Note on storage: *Store all components except the microplate and assay buffer (room temperature) at -70°C for the highest stability. The MMP-3 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-3 enzyme is stable for 4 freeze/thaw cycles. To minimize the number of freeze/thaw cycles, aliquot the MMP-3 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-P126-9090, BML-P127-9090, and BML-PI115-9090 to RT to thaw DMSO.
2. Dilute inhibitor (NNGH, BML-PI115-9090) 1/200 in assay buffer BML-KI127 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. Thaw the DMSO stock vial of substrate BML-P126-9090 and dilute sufficient volume to 40 μ M in assay buffer (10 μ l needed per well). Warm to reaction temperature (e.g. 37°C).
4. Dilute MMP-3 enzyme 1/100 in assay buffer to required total volume (20 μ l 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:
 - Calibration = 80 μ l in 3 wells (see step 11)
 - Control (no inhibitor) = 70 μ l
 - Inhibitor NNGH = 50 μ l
 - Test inhibitor = varies (see Table 1, below)

Example of plate:	<u>well#</u>	<u>sample</u>
	A1	Calibration
	B1	Calibration
	C1	Calibration
	D1	Control
	E1	Control
	F1	Inhibitor NNGH
	G1	Inhibitor NNGH
	H1	Test inhibitor
	I1...	Test inhibitor...

6. Allow microplate to equilibrate to assay temperature (e.g. 37°C).
7. Add 20 μ l MMP-3 (diluted in step 4) to the control, inhibitor NNGH, and test inhibitor wells. Final amount of MMP-3 will be 2 U per well (0.02U/ μ l). Remember to **not** add MMP-3 to the calibration wells!
8. Add 20 μ l NNGH inhibitor (diluted in step 2) to the inhibitor NNGH wells only. Final inhibitor concentration = 1.3 μ M.
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. In the meantime, calibrate the fluorescent microplate reader, using Ex/Em=328/420: Prewarm assay buffer to reaction temperature in 3 wells in the microplate, then to each add 10µl BML-P126-9090 substrate peptide to give the concentration to be used in the assay (e.g., for 4 µM final add 10µl 40µM) and mix. When the fluorescent signal is constant, use this reading as the zero (Blank) value in arbitrary fluorescence units (RFUs). Using the same wells, with their mixtures of substrate peptide and buffer, add 10µl calibration standard peptide BML-P127-9090 to give 3 different final molar concentrations ranging between 2 and 10% of the substrate peptide molar concentration (e.g., 80, 200, and 400 nM) and measure their fluorescence. Use these values to build a standard curve relating micromolar BML-P127-9090 concentration (x axis) to RFUs (y axis). The slope of the line is the conversion factor (CF). If multiple concentrations of substrate peptide are used, such as in kinetic determinations, step 11 must be performed for each concentration, due to absorptive quenching by the substrate peptide. Note: this calibration can be done at any time.
12. Start reactions by the addition of 10 µl BML-P126-9090 substrate (diluted and equilibrated to reaction temperature in step 3). Final substrate concentration = 4 µM.
13. Continuously read plates in the fluorescent microplate reader, using Ex/Em=328/420. For example, record data at 1 minute time intervals for 10 minutes.
14. Perform data analysis (see below).

NOTE: Retain microplate for future use of unused wells.

TABLE 1. Example of Samples.

Sample	Assay buffer	MMP-3 (0.1U/µl)	Inhibitor (6.5 µM)	Substrate (40µM)	Total Volume
Control	70 µl	20 µl	0	10 µl	100 µl
Inhibitor NNGH	50 µl	20 µl	20 µl	10 µl	100 µl
Test inhibitor*	X µl	20 µl	Y µl	10 µl	100 µl

*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 µl.

DATA ANALYSIS

Plotting

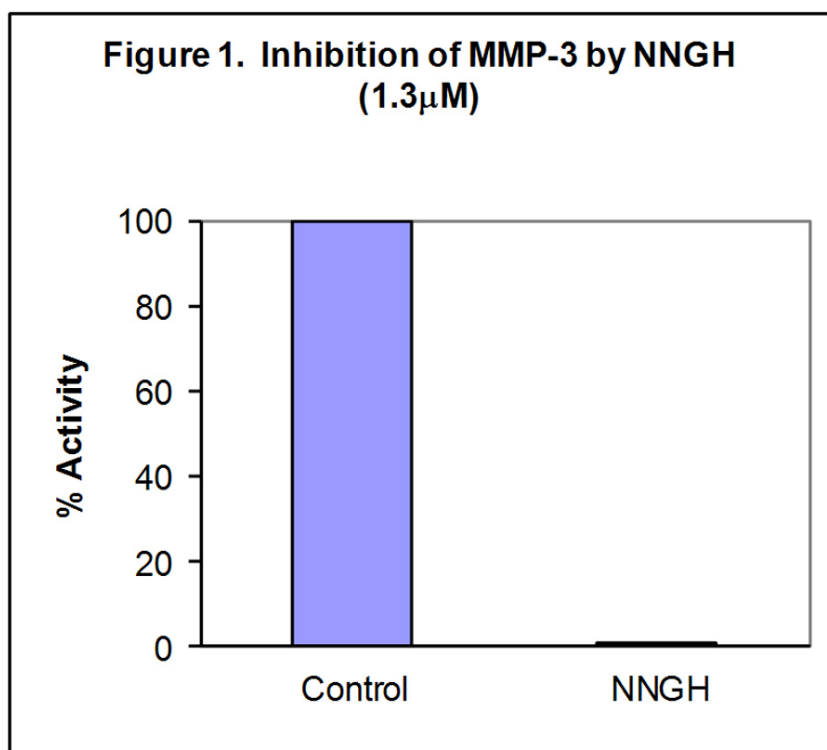
15. Plot data as RFUs (minus Blank RFU value determined during calibration, step 11) versus time for each sample.
16. Determine the range of initial time points during which the reaction is linear.
17. Obtain the initial reaction velocity (**V**) in RFUs/min: determine the slope of a line fit to the initial linear portion of the data plot using an appropriate routine.
18. It is best to use a range of inhibitor concentrations, in duplicate. Average the slopes of duplicate samples.

Data Reduction

To determine inhibitor % remaining activity:

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

See Figure 1 for example of results.



To determine the activity of the samples expressed as picomoles substrate hydrolyzed per minute:

$$X \text{ pmoles substrate/min} = 1/CF \times V \times \text{vol}$$

Where CF is the conversion factor (micromolar concentration/RFUs, from step 11), **V** is initial reaction velocity (RFUs/min, from step 17), and vol is the reaction volume in microliters (100).

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 NONINFRINGEMENT 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.

© 2011



www.enzolifesciences.com
Enabling Discovery in Life

Global Headquarters

Enzo Life Sciences Inc.

10 Executive Blvd

Farmingdale, NY 11735

(p) 1-800-942-0430

(f) 1-631-694-7501

(e) info-usa@enzolifesciences.com

ch@enzolifesciences.com

Enzo Life Sciences (ELS) AG

Industriestrasse 17, Postfach

CH-4415 Lause / Switzerland

(p) +41/0 61 926 89 89

(f) +41/0 61 926 89 79

(e) [info-](mailto:info@enzolifesciences.com)

Please visit our website at www.enzolifesciences.com for additional contact information.