



Enabling Discovery in Life Science®

MMP-12 Fluorometric Drug Discovery Kit, GREEN

Designed to screen MMP-12 inhibitors using a quenched fluorogenic peptide.

**Instruction Manual
BML-AK312**

For research use only

BACKGROUND

Matrix metalloproteinase-12 (MMP-12, metalloelastase, macrophage elastase, commonly confused with neutrophil elastase) 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 specificities¹⁻³. Targets of MMP-12 include elastin, fibronectin, laminin, plasminogen, u-PAR, and tissue factor pathway inhibitor¹⁻⁶. MMP-12 is secreted as a 53 kDa proenzyme (as measured by SDS-PAGE), and activated by cleavage to forms of 45-22 kDa⁷. MMP-12 is an important target for inhibitor screening due to its involvement in diseases such as cancer, allergy and emphysema^{1-3, 8}.

The *MMP-12 Fluorescent Drug Discovery Kit, GREEN* is a complete assay system designed to screen MMP-12 inhibitors using a quenched MMP-3 Fluorogenic Substrate: 5-FAM-Arg-Pro-Lys-Pro-Val-Glu-Nva-Trp-Arg-Lys(TQ2W)-NH₂ [5-FAM = 5-carboxyfluorescein; Nva=norvaline; TQ2W=quencher]. FAM fluorescence is thoroughly quenched by the TQ2W group until cleavage by MMPs separates the two moieties.

The green Fluorogenic MMP substrate provided in the kit, which is cleaved by both MMP-3 and MMP-12, offers key advantages over other MMP substrates. 1) Emission in the green region of the spectrum avoids the interference at lower wavelengths often exhibited by screening compounds, and by substances commonly found in biological samples and tissue culture medium. 2) The ultra-strong fluorescence of this substrate allows for substrate concentrations much lower than the K_m , a condition generally desirable in inhibitor screening/kinetics assays.

The assays are performed in a convenient 96-well microplate format. The kit is useful to screen inhibitors of MMP-12, a potential therapeutic target. The compound NNGH⁷ is also included as a prototypic control inhibitor.

Please contact Enzo Life Sciences for kit components in bulk.

REFERENCES:

1. *Matrix metalloproteinases: they're not just for matrix anymore!*: L.J. McCawley and L.M. Matrisian; *Curr. Opin. Cell Biol.* **13**, 534 (2001)
2. *Updated biological roles for matrix metalloproteinases and new "intracellular" substrates revealed by degradomics*: G.S. Butler and C.M. Overall; *Biochemistry* **48**, 10830 (2009)
3. *Matrix metalloproteinases: regulators of the tumor microenvironment*: K. Kessenbrock and Z. Werb; *Cell* **141**, 52 (2010)
4. *Matrix metalloproteinase degradation of extracellular matrix: biological consequences*: S.D. Shapiro; *Curr. Opin. Cell Biol.* **10**, 602 (1998)
5. *Proteolysis of the urokinase-type plasminogen activator receptor by metalloproteinase-12: implication for angiogenesis in fibrin matrices*. P. Koolwijk *et al.* *Blood* **97**, 3123 (2001)
6. *Matrix metalloproteinases cleave tissue factor pathway inhibitor. Effects on coagulation* A. Belaaouaj *et al.* *J. Biol. Chem.* **275**, 27123 (2000)
7. *Hydrolysis of a broad spectrum of extracellular matrix proteins by human macrophage elastase*. T.J. Gronski, Jr. *et al.* *J. Biol. Chem.* 1997 **272**, 12189 (1997)
8. *Matrix metalloproteinase-12 is a therapeutic target for asthma in children and young adults*. Mukhopadhyay S. *et al.*; *J Allergy Clin Immunol* **126**, 70 (2010).

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.

****DO NOT PERFORM THIS ASSAY IN WHITE PLATE DUE TO PLATE AUTOFLUORESCENCE AT THE WAVELENGTHS USED****

COMPONENTS OF BML-AK312 KIT

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

FORM: *E. coli* recombinant human MMP-12 catalytic domain (calculated MW 20.3 kDa). Purity >95% by SDS-PAGE.

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: 140 U

PRESENTATION: Frozen liquid in screw-cap microfuge vial.

BML-P278-9090 SUBSTRATE (MMP-3 Fluorogenic Substrate; peptide MW=2156.9)

FORM: 250 μ M in DMSO (dimethylsulfoxide)

STORAGE: -70°C

PRESENTATION: 40 μ l in amber screw-cap microfuge vial.

BML-KI583-0010 5-FAM CALIBRATION STANDARD

FORM: 10 μ M in DMSO; MW=376.3

STORAGE: -70°C

PRESENTATION: 20 μ l in amber screw-cap microfuge vial.

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

FORM: 1.3mM in DMSO

STORAGE: -20°C

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

BML-KI175-0020 ASSAY BUFFER

50mM HEPES, 10mM CaCl₂, 0.05% Brij-35, pH 7.5

FORM: Liquid in screw-cap plastic bottle

STORAGE: Room temperature

QUANTITY: 20ml

80-2409 96-WELL BLACK NBS MICROPLATE

STORAGE: Room temperature.

OTHER MATERIALS REQUIRED

1. Fluorescent microplate reader capable of excitation at 494 nm and emission at 521 nm.
2. Pipetmen or multi-channel pipetmen capable of pipetting 1-100 μ l accurately
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 and assay buffer (room temperature) at -70°C for the highest stability. The MMP-12 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-12 enzyme is stable for 4 freeze/thaw cycles. To minimize the number of freeze/thaw cycles, aliquot the MMP-12 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-P278-9090, BML-KI583-0010, and BML-PI115-9090 to RT to thaw DMSO.
2. Dilute inhibitor (NNGH, BML-PI115-9090) 1/200 in assay buffer BML-KI583 as follows. Add $1\mu\text{l}$ inhibitor into $200\mu\text{l}$ assay buffer, in a separate tube.
3. Thaw the DMSO stock vial of substrate BML-P278-9090 and dilute sufficient volume to $5\mu\text{M}$ in assay buffer ($10\mu\text{l}$ needed per well).
4. Dilute MMP-12 enzyme to $50\text{ mU}/\mu\text{l}$ in assay buffer to required total volume ($20\mu\text{l}$ 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 = $80\mu\text{l}$ in 3 wells (see step 11)
Control (no inhibitor) = $70\mu\text{l}$
Inhibitor NNGH = $50\mu\text{l}$
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
	A2...	Test inhibitor

6. Allow microplate to equilibrate to assay temperature (e.g. 37°C).
7. Add $20\mu\text{l}$ MMP-12 (diluted in step 4) to the control, inhibitor NNGH, and test inhibitor wells. Final amount of MMP-12 will be 1 U per well ($10\text{ mU}/\mu\text{l}$).
8. Add $20\mu\text{l}$ NNGH inhibitor (diluted in step 2) to the inhibitor NNGH wells only. Final inhibitor concentration = $1.3\mu\text{M}$.
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=494/521: Prewarm 80µl assay buffer to reaction temperature in 3 wells in the microplate, then to each add 10µl BML-P278-9090 substrate peptide to give the concentration to be used in the assay (e.g., for 0.5µM final add 10µl 5µM) 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-KI583 to give 3 different final molar concentrations ranging between 2 and 10% of the substrate peptide molar concentration (e.g., 10, 25, and 50 nM) and measure their fluorescence. Use these values to build a standard curve relating micromolar BML-KI583 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 10µl BML-P278-9090 substrate (diluted and equilibrated to reaction temperature in step 3). Final substrate concentration = 0.5µM.
13. Continuously read plates in the fluorescent microplate reader, for example, using Ex/Em=494/521 nm, with cutoff ~510 nm. Record data at 1 minute time intervals for 10 minutes at set reaction temperature (e.g. 37°C).
14. Perform data analysis (see below).

NOTE: Retain microplate for future use of unused wells.

TABLE 1. Example of Samples.

Sample	Assay buffer	MMP-12 (50 mU/µl)	Inhibitor (6.5µM)	Substrate (5µM)	Total Volume
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.

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.

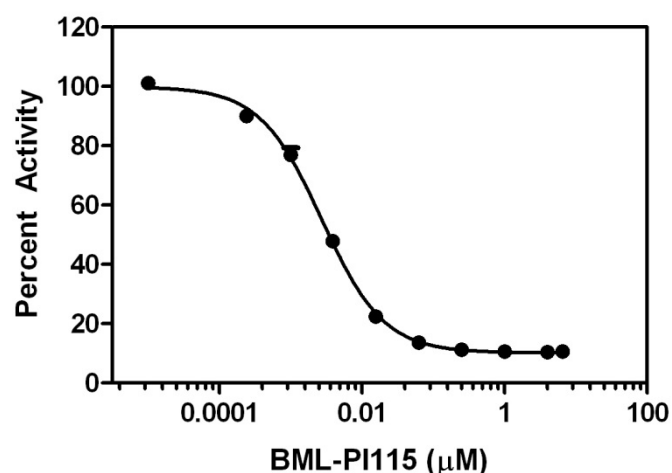


Figure 1. Inhibitor dose response curve: NNGH was pre-incubated with MMP-12 enzyme before reactions were started by the addition of substrate. Final concentrations of reagents were 10 mU/µl MMP-12, 750 nM MMP-3 Fluorogenic Substrate, 0.5 % DMSO. $IC_{50} = 3$ nM. Data were collected on a Biotek Synergy2 plate reader. The dose-response curve was derived from a fit to a four parameter Hill-Slope model, $Y = \text{Bottom} + (\text{Top} - \text{Bottom}) / (1 + 10^{((\text{Log}IC_{50} - X) * \text{HillSlope}))}$ using Graphpad Prism software. Data are presented as mean of duplicate wells \pm SEM.

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)

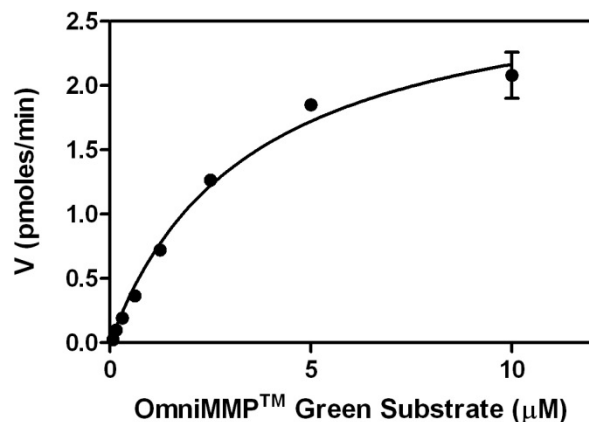


Figure 2. Example graph for k_{cat}/K_m determination for MMP-12. $K_m = 3.5\mu\text{M}$, $k_{cat}/K_m = 1.1 \times 10^5 \text{ M}^{-1}\text{sec}^{-1}$. Data shown are the mean of duplicate wells \pm SEM.

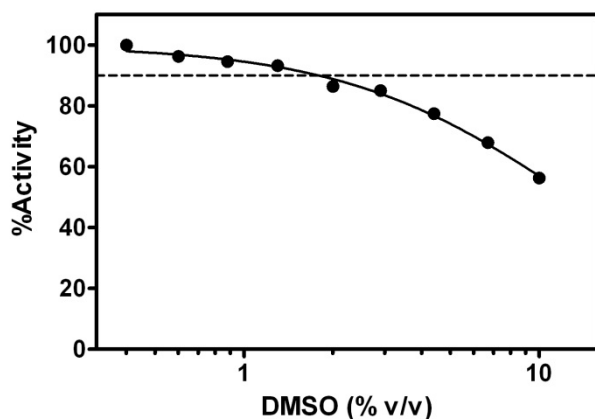


Figure 3. DMSO Sensitivity: Final concentrations of reagents were 10 mU/ μl MMP-12 enzyme, 1 μM MMP-3 Fluorogenic Substrate, variable DMSO concentration. Data are presented as mean of duplicate wells.

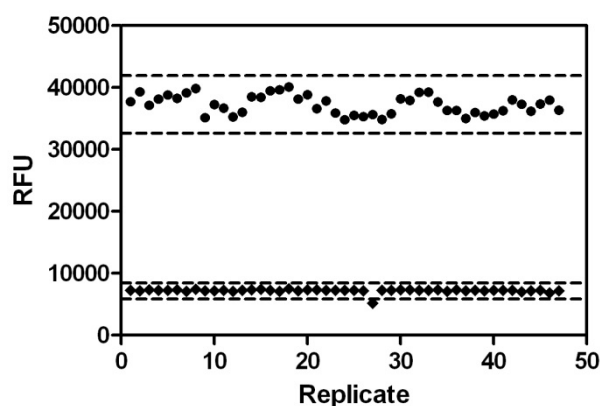


Figure 4. Z analysis: MMP-12 enzyme (●) or BML-KI175 buffer (◆) was prewarmed to 37°C before the addition of MMP-3 Fluorogenic Substrate. After the addition of substrate, the reaction proceeded for 45 minutes before the fluorescence was measured at 485 nm/528 nm (510 nm cutoff) on a BioTek Synergy2 plate reader. Final concentrations were 500 nM substrate, 10 mU/ μl MMP-12, 0.5 % DMSO. The Z factor for this assay was 0.80, ($Z \text{ factor} = 1 - ((3SD^{\text{enzyme}} + 3SD^{\text{buffer}}) / (\text{mean}^{\text{enzyme}} - \text{mean}^{\text{buffer}}))$). Dashed lines indicate 3 * standard deviation.



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.

www.enzolifesciences.com
Enabling Discovery in Life Science[®]

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

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-ch@enzolifesciences.com

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