



# Product Manual

## **MMP-9 Fluorometric Drug Discovery Kit, RED**

Designed to screen MMP-9 inhibitors using a quenched fluorogenic peptide.

**Instruction Manual**

**BML-AK306**

*For research use only*

**✦ MMP-9 Fluorometric Drug Discovery Kit, RED – BML-AK306 ✦****BACKGROUND**

Matrix metalloproteinase-9 (MMP-9, gelatinase B, 92kDa type IV 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>1-3</sup>. Targets of MMP-9 include native and denatured collagens, fibronectin, elastin, laminin, pro-TNF- $\alpha$ , and interleukins and their receptors<sup>1-4</sup>. MMP-9 is secreted as a 92kDa proenzyme (as measured by SDS-PAGE), and activated by cleavage to forms 82kDa and smaller<sup>5</sup>. MMP-9 is an important target for inhibitor screening due to its involvement in diseases such as alopecia<sup>6</sup>, cancer, angiogenesis, and metastasis<sup>3,6,7</sup>.

The *MMP-9 Fluorometric Drug Discovery Kit, RED* is a complete assay system designed to screen MMP-9 inhibitors using a quenched fluorogenic substrate OmniMMP™ RED: TQ3-GABA-Pro-Cha-Abu-Smc-His-Ala-Dab(6-TAMRA)-Ala-Lys-NH<sub>2</sub> [TQ3=quencher; GABA=4-aminobutyric acid; Cha=L-cyclohexylalanine; Abu=2-aminobutyric acid; Smc=S-methyl-L-cysteine; Dab=2,4-diaminobutyric acid; 6-TAMRA=6-tetramethylrhodamine]. TAMRA fluorescence is thoroughly quenched by the TQ3 group until cleavage by MMPs separates the two moieties.

The OmniMMP™ RED substrate offers key advantages over other MMP substrates. 1) Emission at the red end of the spectrum (576nm after excitation at 545nm) avoids the interference at lower wavelengths often exhibited by screening compounds, and by substances commonly found in biological samples and tissue culture medium. 2) MMP substrate peptides display poor aqueous solubility, often with  $K_m$ s near their limits of solubility, making enzyme and inhibitor kinetics difficult. MMP  $K_m$ s for OmniMMP™ RED substrate are well below its solubility limit. 3) OmniMMP™ RED is avidly cleaved by MMPs, with  $k_{cat}/K_m$ s in the range of  $\sim 10^4$ - $10^6$  M<sup>-1</sup>sec<sup>-1</sup>. 4) The ultra-strong fluorescence of OmniMMP™ RED allows for substrate concentrations much lower than the  $K_m$ , a condition generally desirable in inhibitor screening assays.

The assays are performed in a convenient 96-well microplate format. The kit is useful to screen inhibitors of MMP-9, a potential therapeutic target. The compound NNGH<sup>8</sup> 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. *A novel role of metalloproteinase in cancer-mediated immunosuppression*: B.-C. Sheu *et al.*; Cancer Res. **61**, 237 (2001)
5. *Analysis of the role of the COOH-terminal domain in the activation, proteolytic activity, and tissue inhibitor of metalloproteinase interactions of gelatinase B*: J. P. O'Connell *et al.*; J. Biol. Chem. **269**, 14967 (1994)
6. *Identification of clustered cells in human hair follicle responsible for MMP-9 gelatinolytic activity: consequences for the regulation of hair growth*: F. Jarrousse *et al.*; Int. J. Dermatol. **40**, 385 (2001)
7. *Matrix metalloproteinase-9 triggers the angiogenic switch during carcinogenesis*: G. Bergers *et al.*; Nat. Cell Biol. **2**, 737 (2000)
8. *Discovery of CGS 27023A, a non-peptidic, potent, and orally active stromelysin inhibitor that blocks cartilage degradation in rabbits*: L.J. MacPherson *et al.*; J. Med. Chem. **40**, 2525 (1997)

**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 A WHITE PLATE DUE TO PLATE AUTOFLUORESCENCE AT THE WAVELENGTHS USED\*\***

### **COMPONENTS OF BML-AK306 KIT**

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

FORM: *E. coli*. Recombinant human MMP-9 catalytic domain (calculated MW 39kDa) 2.68U/μl

UNIT DEFINITION: One unit is defined as the amount of enzyme that will hydrolyze 100μM thiopeptide Ac-PLG-[2-mercapto-4-methyl-pentanoyl]-LG-OC<sub>2</sub>H<sub>5</sub> (Cat. # BML-P125) at 100 pmol/min @ 37°C.

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

QUANTITY: 130 U

#### **BML-P277-9090 SUBSTRATE (OmniMMP™ RED fluorogenic substrate peptide; MW=1910.7)**

FORM: 250μM in DMSO

STORAGE: -70°C

QUANTITY: 40μl

**BML-KI582-0010 6-TAMRA CALIBRATION STANDARD**FORM: 10 $\mu$ M in DMSO; MW=430.4

STORAGE: -70°C

QUANTITY: 20 $\mu$ l**BML-PI115-9090 INHIBITOR (NNGH; MW=316.4)**

FORM: 1.3 mM in DMSO

STORAGE: -70°C

QUANTITY: 50 $\mu$ l**BML-KI175-0020 ASSAY BUFFER**50mM HEPES, 10mM CaCl<sub>2</sub>, 0.05% Brij-35, pH 7.5

FORM: Liquid in screw-cap plastic bottle

STORAGE: -20°C

QUANTITY: 20ml

**80-2409 96-WELL BLACK NBS MICROPLATE**

STORAGE: Room temperature.

**OTHER MATERIALS REQUIRED**

- Fluorescent microplate reader capable of excitation at 545nm and emission at 576nm. The following Ex/Em has also been used: 530/590.
- Pipetmen or multi-channel pipetmen capable of pipetting 1-100 $\mu$ l accurately.
- Ice bucket to keep reagents cold until use.
- Water bath or incubator for component temperature equilibration.

**EXPERIMENTAL METHODS**

Note on storage: Store all components except the microtiter plate at -70°C for the highest stability. Components with storage temperatures other than -70°C can be stored at the temperature listed OR at -70°C without any deleterious effects. The MMP-9 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-9 enzyme is stable for at least 5 freeze/thaw cycles. To minimize the number of freeze/thaw cycles, aliquot the MMP-9 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-P277-9090, BML-KI582, and BML-PI115-9090 to RT to thaw DMSO.
2. Dilute inhibitor (NNGH, BML-PI115-9090) 1/200 in assay buffer BML-KI582 as follows. Add 1 $\mu$ l inhibitor into 200 $\mu$ l assay buffer, in a separate tube. Warm to reaction temperature (e.g. 37°C).
3. Thaw the DMSO stock vial of substrate BML-P277-9090 and dilute sufficient volume to 5 $\mu$ M in assay buffer (10 $\mu$ l needed per well). Warm to reaction temperature (e.g. 37°C).
4. Dilute MMP-9 enzyme to 45mU/ $\mu$ l in assay buffer to required total volume (20 $\mu$ 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 black microplate as follows:

Calibration = 80 $\mu$ l in 3 wells (see step 11)

Control (no inhibitor) = 70 $\mu$ l

Inhibitor NNGH = 50 $\mu$ l

Test inhibitor = varies (see Table 1)

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 $\mu$ l MMP-9 (diluted in step 4) to the control, inhibitor NNGH, and test inhibitor wells. Final amount of MMP-9 will be 0.9U per well (9mU/ $\mu$ l).
8. Add 20 $\mu$ l NNGH inhibitor (diluted in step 2) to the inhibitor NNGH wells only. Final inhibitor concentration = 1.3 $\mu$ M. Note: 1.3 $\mu$ M NNGH will inhibit MMP-9 by > 96% under these conditions (see **Figure 1**).
9. Add desired volume of test inhibitor to appropriate wells. See **Table 1**. Final DMSO concentrations above 3% have been shown to decrease the rate of MMP-9 in this assay (**Figure 3**).

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=545/576nm, with cutoff set at 570nm: Prewarm 80µl assay buffer to reaction temperature in 3 wells in the microplate, then to each add 10µl BML-P277-9090 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 (diluted in assay buffer) calibration standard BML-KI582 to give 3 different final molar concentrations ranging between 2 and 10% of the substrate peptide molar concentration (e.g., 10, 25, and 50nm) and measure their fluorescence. Use these values to build a standard curve relating micromolar BML-KI582 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-P277-9090 substrate (diluted and equilibrated to reaction temperature in step 3). Final substrate concentration = 0.5µM.
13. Continuously read plates in the fluorescent microplate reader, using Ex/Em=545/576nm, with cutoff set at 570nm. For example, 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.

<b>Sample</b>	<b>Assay buffer</b>	<b>MMP-9 (45mU/µl)</b>	<b>Inhibitor (6.5µM)</b>	<b>Substrate (5µM)</b>	<b>Total Volume</b>
<b>Control</b>	70µl	20µl	0	10µl	100µl
<b>Inhibitor NNGH</b>	50µl	20µl	20µl	10µl	100µl
<b>Test inhibitor*</b>	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

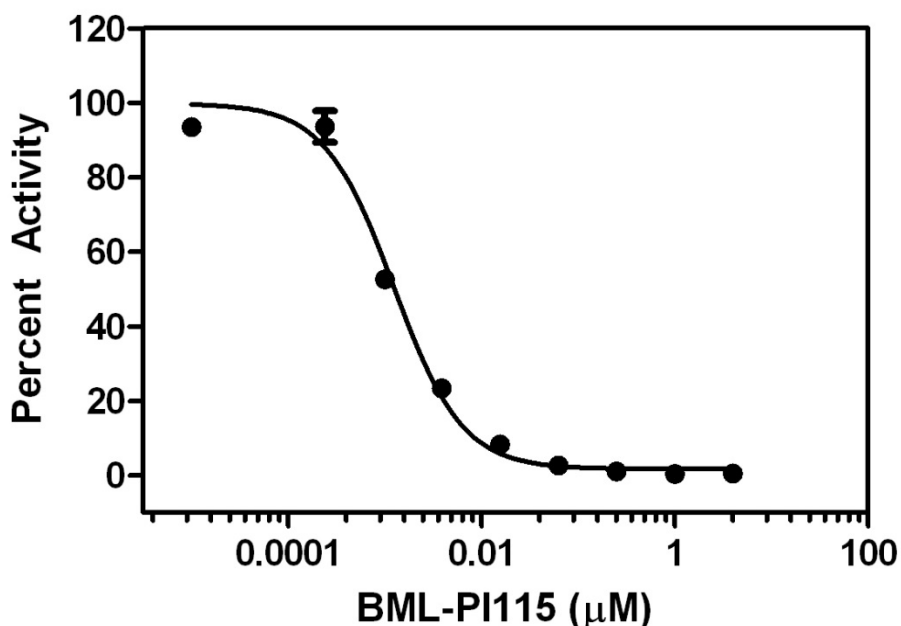
15. Plot data as RFUs (minus Blank RFU value determined during calibration, step 11) versus time for each sample.
16. Determine the range of initial time points during which the reaction is linear.
17. Obtain the initial reaction velocity (**V**) in RFUs/min: determine the slope of a line fit to the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

### Data Reduction

To determine inhibitor % remaining activity:

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

See **Figure 1** for example of results.

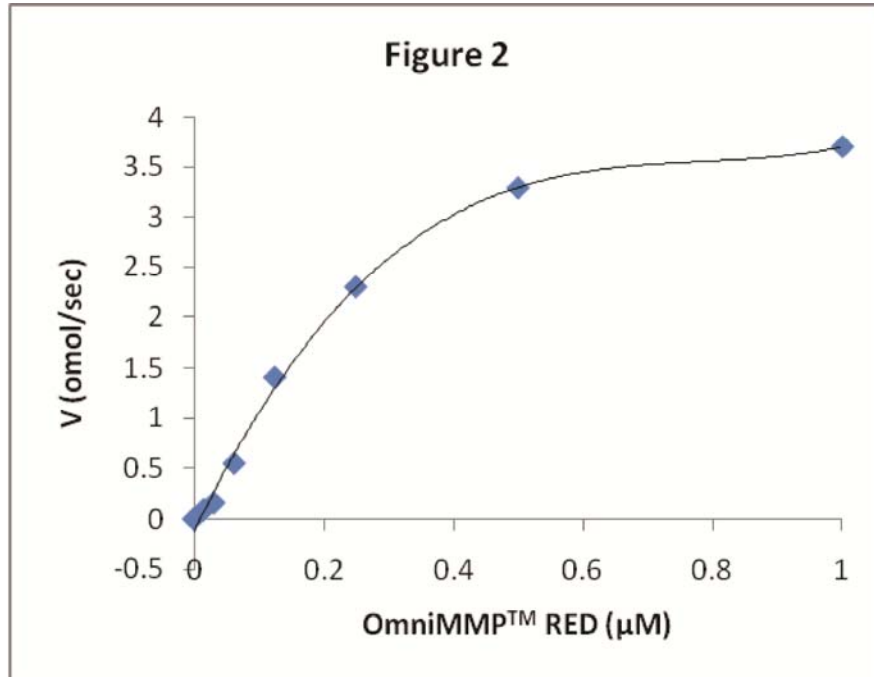


**Figure 1.** Inhibitor dose response curve: NNGH was pre-incubated with MMP-9 enzyme before reactions were started by the addition of substrate. Final concentrations of reagents were 9 mU/µl MMP-9, 500nm OmniMMP™ Red, 1% DMSO. IC<sub>50</sub> was calculated as 1.2nm. Data are presented as mean of duplicate wells ± SEM.

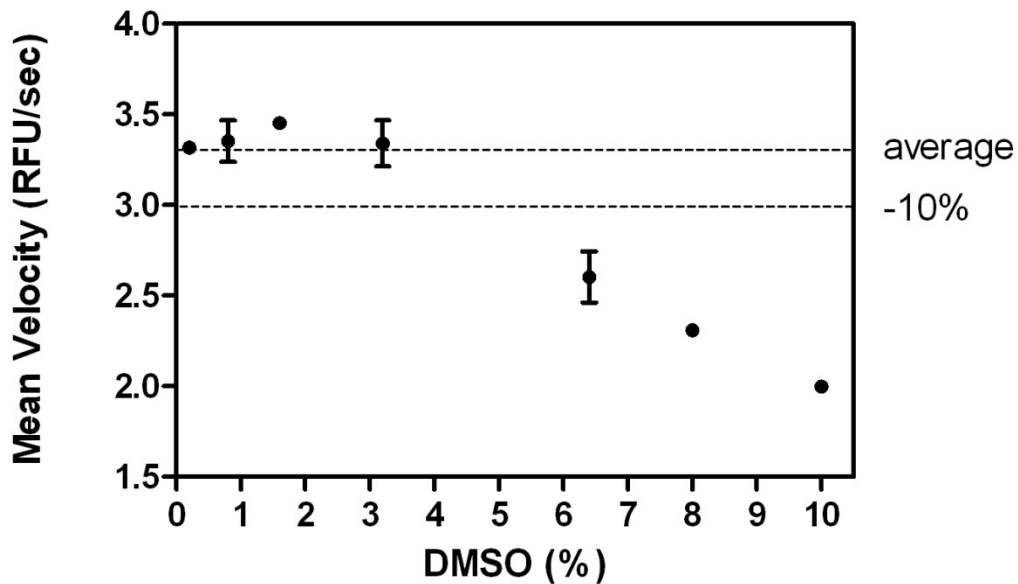
To determine the activity of the samples expressed as picomoles substrate hydrolyzed per minute:

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

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).

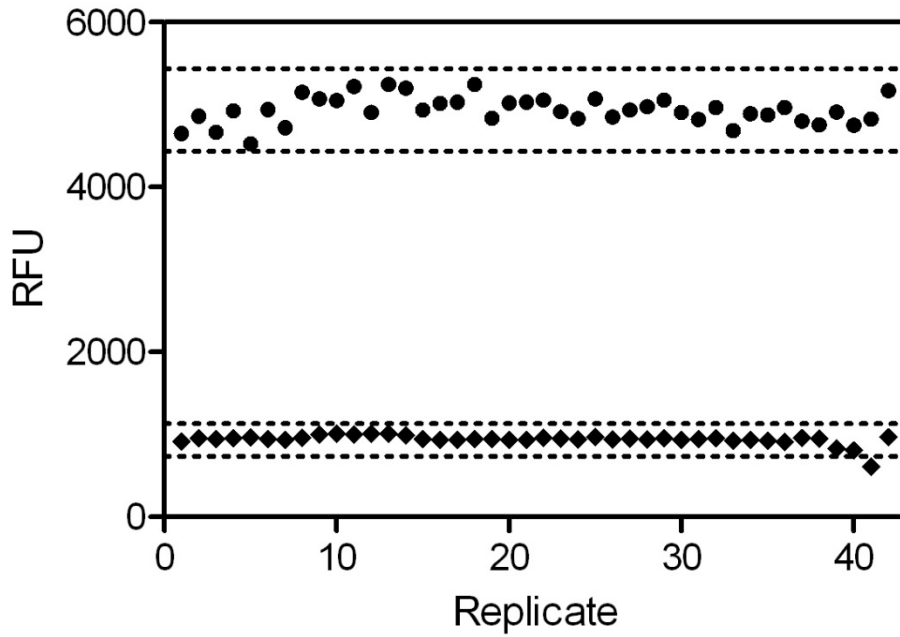


**Figure 2.** Example graph for  $k_{cat}/K_m$  determination for MMP-9.  $K_m = 0.35 \mu\text{M}$ ,  $k_{cat}/K_m = 2.3 \times 10^7 \text{ M}^{-1}\text{sec}^{-1}$ .



**Figure 3.** DMSO Sensitivity: Final concentrations of reagents were 9 mU/μl MMP-9 enzyme, 500nm OmniMMP™ Red, variable DMSO concentration. Data are presented as mean of duplicate wells ± SEM.





**Figure 4.** Z' analysis: MMP-9 enzyme was pre-incubated without (●) or with (◆) NNGH Inhibitor for 30 minutes at ~25°C before the addition of OmniMMP™ Red substrate. After the addition of substrate, the reaction proceeded at ~25°C before the fluorescence was measured at 530nm/590nm, 515nm cutoff on a BioTek Synergy2 plate reader. Final concentrations were 500nm substrate, 9 mU/μl MMP-9, 1μM NNGH, 0.5% DMSO. The Z' factor for this assay was 0.83, ( $Z' \text{ factor} = 1 - ((3SD^{\text{control}} + 3SD^{\text{inhibitor}}) / (\text{mean}^{\text{control}} - \text{mean}^{\text{inhibitor}}))$ ). Dashed lines indicate 3 \* standard deviation.



# Product Manual

Notes



## **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.



# 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](http://www.enzolifesciences.com)