

# MMP-3 fluorogenic substrate

Long emission wavelength reduces background

Highly-quenched, ultra-bright fluorogenic substrate for MMP-3 (other MMPs not tested, but likely). 5'-FAM fluorescence is thoroughly quenched by the TQ2W group until cleavage by MMP-3 separates the two moieties.

## Ordering Information

[Order Online »](#)

BML-P278-0100	0.1mg
---------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Avoid freeze/thaw cycles of solution. Protect from light and moisture.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Appearance	Lyophilized orange or red solid.
Application Notes	Useful for inhibitor screening, kinetic analysis, and cellular activity assay.
MW	2156.9
Purity	≥90% (HPLC)
Sequence	5'-FAM-Arg-Pro-Lys-Pro-Val-Glu-Nva-Trp-Arg-Lys(TQ2W)-NH <sub>2</sub>
Solubility	Soluble in DMSO (5mM).
Technical Info / Product Notes	<p>This substrate offers key advantages over other MMP-3 substrates.</p> <ol style="list-style-type: none"><li>1. Emission at the green end 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.</li><li>2. MMP substrate peptides display poor aqueous solubility, often with <math>K_m</math>s near or above their limits of solubility, making enzyme and inhibitor kinetics difficult. MMP-3 <math>K_m</math> (~0.64μM) for this substrate is below its solubility limit (~2μM in assay buffer), allowing for substrate concentrations higher than the <math>K_m</math>, a condition generally desirable in endpoint assays.</li><li>3. In addition to the efficient binding as exhibited by low <math>K_m</math>, this substrate is avidly cleaved by MMP-3, with a <math>k_{cat}/K_m</math> of <math>\sim 6.5 \times 10^7 M^{-1} sec^{-1}</math>.</li><li>4. The high <math>k_{cat}/K_m</math> and the ultra-strong fluorescence of this substrate allow for substrate concentrations much lower than the <math>K_m</math>, a condition generally desirable in inhibitor screening/kinetics assays.</li></ol>



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)