

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 ₂
Solubility	Soluble in DMSO (5mM).

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.
2. MMP substrate peptides display poor aqueous solubility, often with K_m s near or above their limits of solubility, making enzyme and inhibitor kinetics difficult. MMP-3 K_m (~0.64 μ M) for this substrate is below its solubility limit (~2 μ M in assay buffer), allowing for substrate concentrations higher than the K_m , a condition generally desirable in endpoint assays.
3. In addition to the efficient binding as exhibited by low K_m , this substrate is avidly cleaved by MMP-3, with a k_{cat}/K_m of $\sim 6.5 \times 10^7 \text{M}^{-1}\text{sec}^{-1}$.
4. The high k_{cat}/K_m and the ultra-strong fluorescence of this substrate allow for substrate concentrations much lower than the K_m , a condition generally desirable in inhibitor screening/kinetics assays.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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