

TACE substrate (fluorogenic)

Fluorogenic substrate for TACE (ADAM17) (between Ala-Val), as well as MMP-3, MMP-7, MMP-12, and MMP-17, but not for ADAM19, MMP-13, nor MMP-14. Similar substrates are also cleaved by ADAM10, ADAM8 and ADAM9, and MMP-1, -2, and -9. This quenched substrate is useful for TACE inhibitor screening and kinetic analysis. Ex: 328 nm, Em: 393 nm, although the following Ex/Em have been used: 320-340/400-420. Also available: OMNIMMP® Fluorogenic Control peptide (BML-P127).

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P132-0500	0.5mg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name ADAM17 substrate

CAS 192723-42-5

Formula $C_{69}H_{103}N_{23}O_{24}$

MW 1638.7

Purity ≥96%

Sequence Mca-Pro-Leu-Ala-Gln-Ala-Val-Dpa-Arg-Ser-Ser-Ser-Arg-NH₂
[Mca=(7-methoxycoumarin-4-yl)acetyl; Dpa=N-3(2,4-dinitrophenyl)-L-α,β-diaminopropionyl]

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