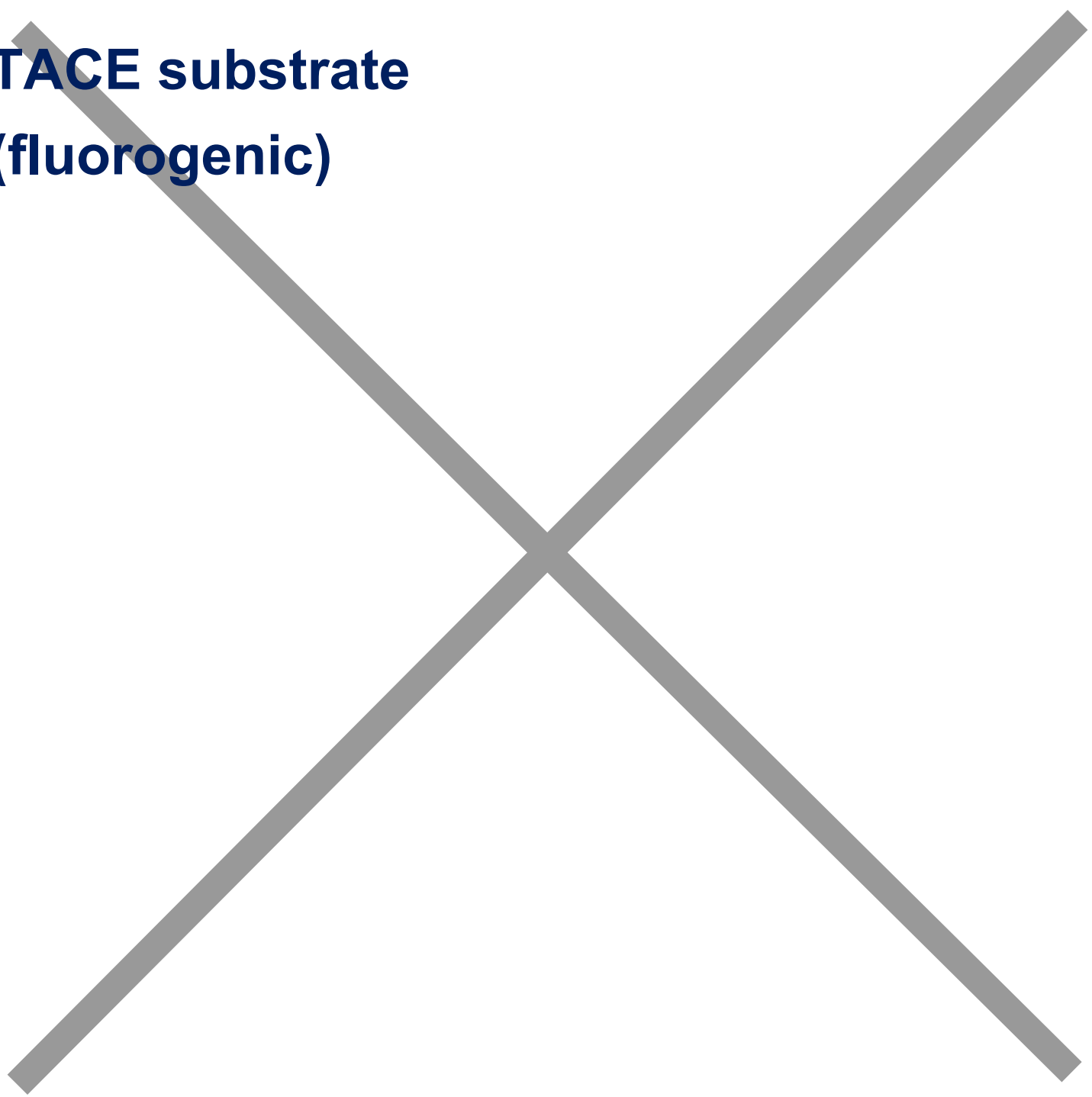


TACE substrate (fluorogenic)



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-NH ₂ [Mca=(7-methoxycoumarin-4-yl)acetyl; Dpa=N-3(2,4-dinitrophenyl)-L-α,β-diaminopropionyl]

Last modified: May 29, 2024