

TACE substrate II (fluorogenic)

Sensitive fluorogenic substrate for TACE (TNF α converting enzyme, ADAM17). Also cleaved by ADAM10 and likely other ADAMs and MMPs. The k_{cat}/K_m of TACE for this substrate is 7-fold higher than for the same peptide with the Mca-Dpa fluorophore-quencher pair. FAM fluorescence is quenched by the TAMRA group until cleavage separates them. Ex.: 485nm, Em.: 535nm, although the following Ex/Em can also be used: 490,494/515,520nm. This substrate is useful for inhibitor screening, kinetic analysis, and cellular activity assay.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P228-0100	100 μ g
---------------	-------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	ADAM17 substrate II (fluorogenic)
Appearance	Solid.
Formulation	Lyophilized.
MW	2157.3
Purity	≥95% (HPLC)
Sequence	5-FAM-Ser-Pro-Leu-Ala-Gln-Ala-Val-Arg-Ser-Ser-Ser-Arg-Lys(5-TAMRA)-NH ₂
Solubility	Soluble in DMSO.



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)