

# ADAM substrate (fluorogenic)

Fluorogenic substrate for ADAM8 ( $k_{\text{cat}}/K_{\text{m}} = 1.0 \times 10^5 \text{ M}^{-1}\text{s}^{-1}$ ), ADAM10 ( $6.2 \times 10^3$ ), ADAM12 ( $2.8 \times 10^5$ ), TACE (TNF- $\alpha$  converting enzyme, ADAM17) ( $4.3 \times 10^5$ ), and likely other ADAMs, especially those that cleave TNF- $\alpha$  type sequences. Also efficiently cleaved by MMP-1, MMP-2, MMP-3, MMP-8, MMP-9, MMP-13, and MMP-14. FAM fluorescence is quenched by the Dabcyl group until cleavage separates them. Ex.: 485nm, Em.: 530nm. This substrate is useful for inhibitor screening and kinetic analysis.

Citations: 6

[View Online »](#)

## Ordering Information

[Order Online »](#)

|               |     |
|---------------|-----|
| BML-P235-0001 | 1mg |
|---------------|-----|

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

|                   |                                                                                                                            |
|-------------------|----------------------------------------------------------------------------------------------------------------------------|
| Use/Stability     | As indicated on product label or CoA when stored as recommended. Stock solutions in DMSO are stable for 3 months at -20°C. |
| Handling          | Protect from light and moisture.                                                                                           |
| Long Term Storage | -20°C                                                                                                                      |
| Shipping          | Blue Ice                                                                                                                   |

## Regulatory Status

RUO - Research Use Only

## Product Details

|            |                                                                                                                                              |
|------------|----------------------------------------------------------------------------------------------------------------------------------------------|
| Appearance | Pink powder.                                                                                                                                 |
| MW         | 1542.5                                                                                                                                       |
| Purity     | ≥95% (HPLC)                                                                                                                                  |
| Quantity   | 1mg net peptide.                                                                                                                             |
| Sequence   | Dabcyl-Leu-Ala-Gln-Ala-Homophe-Arg-Ser-Lys(5-FAM)-NH <sub>2</sub> . TFA (Dabcyl=4-(4-dimethylaminophenylazo)benzoyl; FAM=carboxyfluorescein) |
| Solubility | Soluble in DMSO (≥10mM) or TRIS, pH 8.0 assay buffer (≥10μM).                                                                                |

Last modified: May 29, 2024



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)