

# ADAM17 fluorometric drug discovery kit

The *ADAM17 Fluorometric (also known as fluorimetric) Drug Discovery Kit* is a complete assay system designed to screen ADAM17 (TACE) inhibitors using a quenched fluorogenic peptide: Mca-PLAQAV-Dpa-RSSSR-NH<sub>2</sub>. Mca fluorescence is quenched by the Dpa group until cleavage by proteases separates the two moieties. The assays are performed in a convenient 96-well microplate format. The kit is useful to screen inhibitors of ADAM17, a potential therapeutic target. The compound GM6001 is also included as a prototypic control inhibitor.

TACE/ADAM17 (Tumor necrosis factor- $\alpha$ -converting enzyme; A Disintegrin And Metalloproteinase 17) is a soluble or membrane-bound metalloproteinase primarily responsible for activation of proTNF- $\alpha$ , while also targeting proteins such as fractalkine, amyloid precursor proteins, and CD40. ADAM17/TACE is involved in cancer, vascular disorders, and inflammatory diseases such as rheumatoid arthritis and focal ischemic injury. The catalytic domain of ADAM17/TACE is able to cleave proTNF- $\alpha$  and is used in inhibitor screening.

## Ordering Information

[Order Online »](#)

BML-AK310-0001	96 wells
----------------	----------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -80°C

**Shipping** Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** TACE, A disintegrin and metalloproteinase 17

**Application** Activity assay, Fluorescent detection, HTS

**Application Notes** Designed to screen ADAM17 (TACE) inhibitors using a quenched fluorogenic peptide.

**Contents**

- 1 vial ADAM17 enzyme
- 1 vial Substrate (Mca-PLAQAV-Dpa-RSSSR-NH<sub>2</sub>)
- 1 vial CALIBRATION STANDARD
- 1 vial control inhibitor (GM6001)
- 1 bottle (20 ml) assay buffer
- 1 black 96-well microplate
- Instructions

**UniProt ID** P78536



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)