

# ADAM17 (catalytic domain) (human), (recombinant) (His-tag)

ADAM17/TACE 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 can be used in inhibitor screening.

Citations: 5

[View Online »](#)

## Ordering Information

[Order Online »](#)

BML-SE268-0010	10 $\mu$ g
----------------	------------

## Manuals, SDS & CofA

[View Online »](#)

**kDa**

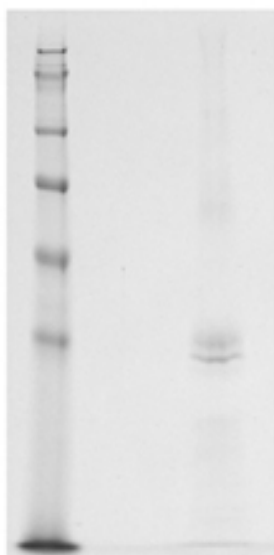
148

98

64

50

36



1

2

SDS-PAGE analysis: Lane 1: MW Marker; Lane 2: 1.0  $\mu$ g of Prod. No. BML-SE268 ADAM17 (catalytic domain) (human), (recombinant) (His-tag).

## Handling & Storage

<b>Use/Stability</b>	Salts (sodium chloride, calcium chloride, etc.) in the assay are inhibitory. ADAM17/TACE is stable after 6 freeze-thaws at ~0.4µg/µl; freeze-thaw stability of more dilute preparations has not been tested and could lead to loss of activity.
<b>Handling</b>	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	TACE, A disintegrin and metalloproteinase 17, Tumor necrosis factor-α-converting enzyme
<b>Application Notes</b>	Useful tool to study enzyme kinetics, cleave target substrates, screen inhibitors.
<b>Formulation</b>	Liquid. In 22.5mM TRIS, pH 7.5, containing 4.5µM ZnCl <sub>2</sub> , 0.0045% Brij-35 and 10% glycerol.
<b>MW</b>	~30.5kDa (calculated), ~36kDa doublet (SDS-PAGE)
<b>Purity</b>	≥90% (SDS-PAGE)
<b>Purity Detail</b>	Purified by multi-step chromatography.
<b>Sequence</b>	Recombinant glycosylated catalytic domain (aa Pro <sup>18</sup> -Val <sup>477</sup> ) of ADAM17/TACE (A disintegrin and metalloproteinase 17; Tumor necrosis factor-α-converting enzyme), cloned from human cDNA (NM_003183), secreted as mature, active enzyme from insect cells, and purified using a C-terminal His-tag.
<b>Source</b>	Produced in insect cells. Produced in a baculovirus expression system.
<b>Specific Activity</b>	≥1800 U/µg enzyme. One unit will hydrolyze one pmole Mca-PLAQAV-Dpa-RSSSR-NH <sub>2</sub> substrate (Prod. No. BML-P132) (10µM) per minute at 37°C, in 25mM TRIS, pH 9.0.

### UniProt ID



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 <a href="mailto:info-usa@enzolifesciences.com">info-usa@enzolifesciences.com</a>	<b>P78536</b> European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 <a href="mailto:info-eu@enzolifesciences.com">info-eu@enzolifesciences.com</a>	Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 <a href="mailto:info-be@enzolifesciences.com">info-be@enzolifesciences.com</a>	France Phone: +33 472 440 655 <a href="mailto:info-fr@enzolifesciences.com">info-fr@enzolifesciences.com</a>	Germany Phone: +49 7621 5500 526 <a href="mailto:info-de@enzolifesciences.com">info-de@enzolifesciences.com</a>	UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 <a href="mailto:info-uk@enzolifesciences.com">info-uk@enzolifesciences.com</a>
---	---	--	--	---	---