## ADAM17 (catalytic domain) (human), (recombinant) (Histag)

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** 

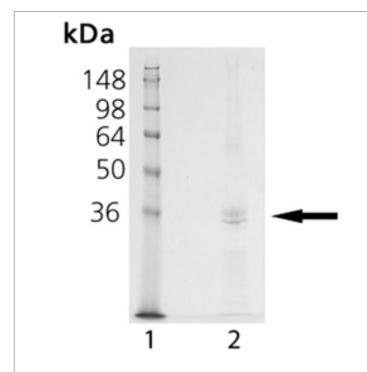
**Order Online** »

BML-SE268-0010

10µg

Manuals, SDS & CofA

**View Online** »



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

## **Handling & Storage**

**Use/Stability** 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.

**Handling** Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.

Long Term Storage -80°C

Shipping Dry Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

Alternative Name TACE, A disintegrin and metalloproteinase 17, Tumor

necrosis factor-α-converting enzyme

Application Notes Useful tool to study enzyme kinetics, cleave target

substrates, screen inhibitors.

Formulation Liquid. In 22.5mM TRIS, pH 7.5, containing 4.5µM ZnCl<sub>2</sub>,

0.0045% Brij-35 and 10% glycerol.

**MW** ~30.5kDa (calculated), ~36kDa doublet (SDS-PAGE)

Purity ≥90% (SDS-PAGE)

Purity Detail Purified by multi-step chromatography.

Sequence Recombinant glycosylated catalytic domain (aa Pro<sup>18</sup>-Val

477) 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.

Source Produced in insect cells, Produced in a baculovirus

expression system.

Specific Activity ≥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 P78536

Last modified: May 29, 2024

