

ADAM17 (CT)

polyclonal antibody

TNF-alpha converting enzyme (TACE) is a membrane-bound metalloprotease-disintegrin in the family of mammalian ADAM (for a disintegrin and metalloprotease), and is responsible for the cleavage of membrane-bound precursor TNF-alpha to its mature, secreted form. TACE also processes other cell surface proteins, including TNF receptor, TGFalpha, the L-selectin adhesion molecule, and alpha-cleavage of amyloid protein precursor (APP).

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 6

[View Online »](#)

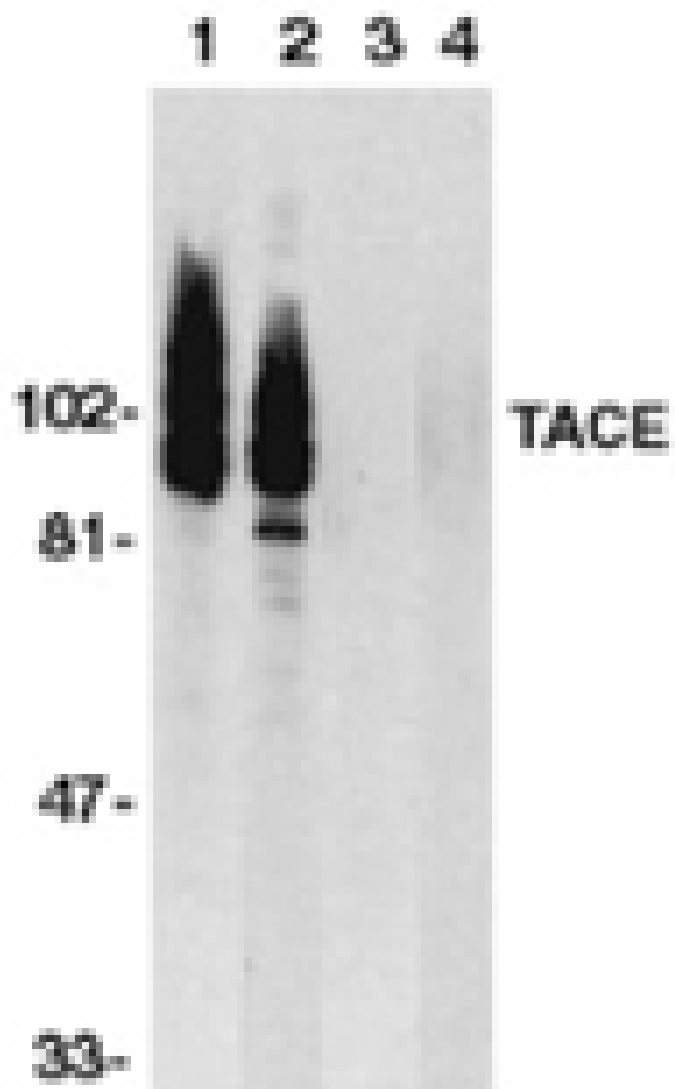
Ordering Information

[Order Online »](#)

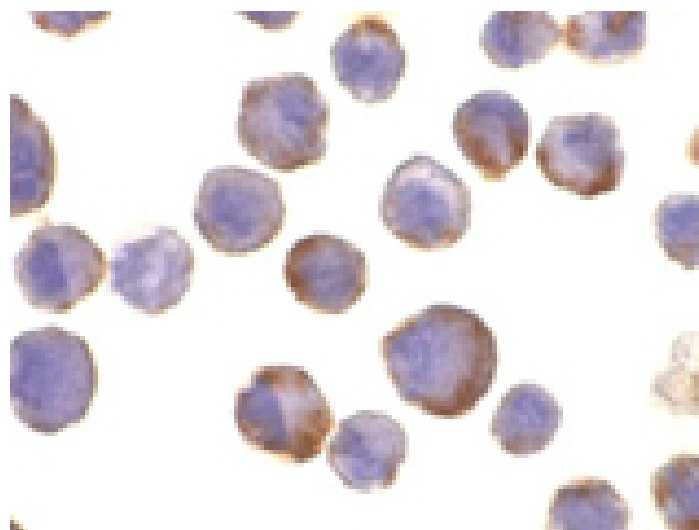
ADI-905-249-100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of TACE in HeLa and Jurkat whole cell lysate in the absence or presence of blocking peptide with TACE antibody at 1:500 dilution.



Immunocytochemistry analysis of TACE in HeLa cells with TACE antibody at 10 μ g/ml.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name TACE, A disintegrin and metalloproteinase 17

Application ICC, WB

Application Notes Detects bands of ~80 to ~130kDa (may represent mature protein, precursor, and glycosylated TACE) by Western blot.

Formulation Liquid. In PBS containing 0.02% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to the sequence near the C-terminus of human ADAM17.

Purity Detail Peptide affinity purified.

Recommendation Dilutions/Conditions Western Blot (1:500-1:2,000) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity Human, Mouse, Rat

UniProt ID P78536

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



Enzo Life Sciences, INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com