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: 5

View Online »

Ordering Information

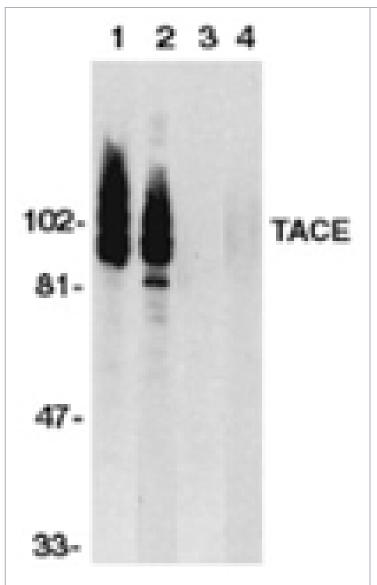
Order Online »

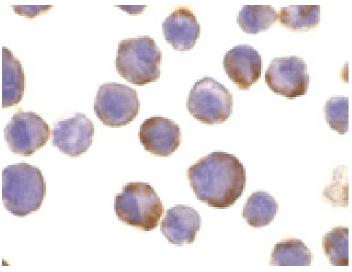
ADI-905-249-100

100µg

Manuals, SDS & CofA

View Online »





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

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.

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 NotesDetects 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/ConditionsWestern 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

•

