

TWEAK antibody (pAb)

Rabbit Anti-Human/Mouse/Rat TNF-like Weak Inducer of Apoptosis (TNF-alpha Related Protein)

Instruction Manual

Catalog Number	PK-AB718-3791
Synonyms	TWEAK Antibody: TNF-related weak inducer of apoptosis, TNF ligand superfamily member 12
Description	TNF-related weak inducer of apoptosis (TWEAK) is a member of the tumor necrosis factor superfamily (TNFSF) of structurally related cytokines. Like most other members of this family, TWEAK is a cell surface-associated type II transmembrane protein although a smaller, biologically active form can also be generated by cleavage near the cell membrane. TWEAK has multiple biological activities, including stimulation of cell growth and angiogenesis, induction of inflammatory cytokines, in addition to stimulation of apoptosis. The TWEAK signal transduction pathway has not been well established but it appears to signal via TweakR/Fn14 in a manner similar to that described for other TNFSF members that bind receptors lacking death domains. Despite its predicted molecular weight of 27kDa, TWEAK may migrate at aberrant positions in SDS-PAGE.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal TWEAK antibody was raised against recombinant human TWEAK (Genbank accession No. NP_003800).
Purification Method	Purified IgG
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	
Formulation	Antibody is supplied in PBS containing 0.02% sodium azide.
Reconstitution	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
Storage & Stability	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB, IHC INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. TWEAK antibody can be used for the detection of TWEAK by Western blot at 1 µg/mL. Antibody can also be used for immunohistochemistry starting at 10 µg/mL.
Images	Available upon request.
References	Yoshida H, Matsui T, Yamamoto T, et al. XBP1 mRNA is induced by ATF6 and spliced by IRE1p in response to ER stress to produce a highly active transcription factor. <i>Cell</i> 2001; 107:881-91. Chicheportiche Y, Bourdon PR, Xu H, et al. TWEAK, a new secreted ligand in the tumor necrosis factor family that weakly induces apoptosis. <i>J. Biol. Chem.</i> 1997; 272:32401-10. Lynch CN, Wang YC, Lund JK, et al. TWEAK induces angiogenesis and proliferation of endothelial cells. <i>J. Biol. Chem.</i> 1999; 274:8455-9. Chicheportiche Y, Chicheportiche R, Sizing I, et al. Proinflammatory activity of TWEAK on human dermal fibroblasts and synoviocytes: blocking and enhancing effects of anti-TWEAK monoclonal antibodies. <i>Arthr. Res.</i> 2002; 4:126-33. 5. Kaplan MJ, Lewis EE, Shelden EA, et al. The apoptotic ligands TRAIL, TWEAK, and Fas ligand mediate monocyte death induced by autologous lupus T cells. <i>J. Immunol.</i> 2002; 169:6020-9.
Images	Available upon request.
Related Products	TWEAK Peptide, Cat.No. PK-AB718-3791P HT-29 Cell Lysate, Cat. No. PK-AB718-1218

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.