

Rabbit Anti-Human Fn14 (Fibroblast Growth Factor Inducible 14, TWEAK Receptor)

Instruction Manual

Catalog Number	PK-AB718-3951
Synonyms	Fn14 Antibody: Fibroblast growth factor-inducible 14, TNFRSF12, TWEAK-R
Description	TWEAK is a TNF family member produced by IFN- γ -stimulated monocytes and induces multiple cell death pathways including caspase-dependent apoptosis and cathepsin B-dependent necrosis. These pathways are mediated by the binding of TWEAK by fibroblast growth factor 14 (Fn14), a member of the TNF receptor superfamily. Yeast two-hybrid experiments have shown that FN14 will bind to the TNFR-associated factors TRAF1, TRAF2, TRAF3, and TRAF5, suggesting that the signaling pathway induced by TWEAK may be mediated by one or more of these adaptor molecules. Fn14 is induced during liver regeneration and is highly expressed in hepatocellular carcinomas. In addition, TWEAK induces liver progenitor cell proliferation, suggesting Fn14 may play a role in hepatocyte growth control and liver neoplasia.
Quantity	100 μ g
Source / Host	Rabbit
Immunogen	Fn14 antibody was raised against a human IgG1 Fc-tagged Fn14 protein.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human
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 μ l 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, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. Fn14 antibody can be used for the detection of Fn14 by Western blot at 2 - 4 μ g/mL. Antibody can also be used for immunohistochemistry starting at 2.5 μ g/mL. For immunofluorescence start at 20 μ g/mL.
Images	Available upon request.
References	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. Nakayama M, Ishidoh K, Kojima Y, et al. Fibroblast growth factor-inducible 14 mediates multiple pathways of TWEAK-induced cell death. <i>J. Immunol.</i> 2003; 170:341-8. Brown SA, Richards CM, Hanscom HN, et al. The Fn14 cytoplasmic tail binds tumour-necrosis-factor-receptor-associated factors 1, 2, 3, and 5 and mediates nuclear factor-kappaB activation. <i>Biochem. J.</i> 2003; 371:395-403. Feng SL, Guo Y, Factor VM, et al. The Fn14 immediate-early response gene is induced during liver regeneration and highly expressed in both human and murine hepatocellular carcinomas. <i>Am. J. Pathol.</i> 2000; 156:1253-61
Images	Available upon request.
Related Products	HepG2 Lysate, Cat. No. PK-AB718-1211 TRAF3 /1 Antibody, Cat. No. PK-AB718-3599 TWEAK Antibody, Cat. No. PK-AB718-3791

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