

TOCA-1 (NT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat TOCA-1 (NT)

Instruction Manual

Catalog Number	PK-AB718-4375
Synonyms	TOCA-1 Antibody; Transducer of cdc42-dependent actin assembly 1, formin binding protein 1-like, FNBP1L
Description	Actin reorganization is important for the regulation of neuronal morphology. A protein involved in this process, the transducer of cdc42-dependent actin assembly 1 (TOCA-1) protein, a member of the evolutionarily conserved pombe CDC15 homology (PCH) protein family, is an essential component of the Cdc42 pathway. TOCA-1 binds both N-WASP and Cdc42 and is essential for Cdc42- and PIP2-induced actin polymerization, suggesting that TOCA-1 mediates Cdc42-dependent actin nucleation by activating the N-WASP-WIP complex. Decreased expression of TOCA-1 significantly enhances neurite elongation in PC-12 cells; its overexpression in the same cells suppresses neurite elongation. It has been suggested that TOCA-1 negatively regulates axon branching by regulating membrane trafficking by regulating membrane trafficking through the F-BAR/EFC domain. Multiple isoforms of TOCA-1 are known to exist.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	TOCA-1 antibody was raised in rabbits against a 15 amino acid peptide near the amino terminus of the human TOCA-1.
Purification Method	Affinity chromatography purified via peptide column.
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, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. TOCA-1 antibody can be used for detection of TOCA-1 by Western blot at 0.5 µg/mL. Antibody can also be used for immunohistochemistry starting at 2.5 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Ho H-Y H, Rohatgi R, Lebensohn AM, et al. Toca-1 mediates Cdc42-dependent actin nucleation by activating the N-WASP-WIP complex. Cell 2004; 118:203-16. Kakimoto T, Katoh H, and Negishi M. Regulation of neuronal morphology by Toca-1, an F-BAR/EFC protein that induces plasma membrane invagination. J. Biol. Chem. 2006; 281:29042-43.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4375P; TOCA-1 Peptide Cat.No. PK-AB718-1303; Human Brain Tissue Lysate

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