

MANF (NT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat MANF (NT)

Instruction Manual

Catalog Number	PK-AB718-4349
Synonyms	MANF Antibody: Mesencephalic astrocyte-derived neurotrophic factor, arginine-rich mutated in early stage tumor, ARMET, ARP
Description	MANF, also known as ARMET, was initially identified as a protein containing an arginine-rich region that was highly mutated in a variety of tumors. More recently it was identified as a mesencephalic astrocyte-derived neurotrophic factor with selectivity for dopaminergic neurons, similar to glial cell line-derived neurotrophic factor (GDNF) and CDNF. In rat brain slices, MANF enhanced nigral gamma-aminobutyric acid release. Like GDNF and CDNF, MANF has selective neuroprotective activity for dopaminergic neurons suggesting that it may be indicated for the treatment of Parkinson's disease. Expression of MANF has also been shown to be induced during ER stress, suggesting that it may play a role in protein quality control during ER stress. This antibody does not cross-react with CDNF.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	MANF antibody was raised in rabbits against a 12 amino acid peptide from near the amino terminus of human MANF.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	This antibody does not cross-react with CDNF.
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 Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. MANF antibody can be used for detection of MANF by Western blot at 1 - 2 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	Shridhar V, Rivard S, Shridhar R, et al. A gene from human chromosomal band 3p21.1 encodes a highly conserved arginine-rich protein and is mutated in renal cell carcinomas. <i>Oncogene</i> 1996; 12:1931-9. Shridhar R, Shridhar V, Rivard S, et al. Mutations in the arginine-rich protein gene, in lung, breast, and prostate cancers, and in squamous cell carcinoma of the head and neck. <i>Cancer Res.</i> 1996; 56:5576-8. Petrova P, Raibekas A, Pevsner J, et al. MANF: a new mesencephalic, astrocyte-derived neurotrophic factor with selectivity for dopaminergic neurons. <i>J. Mol. Neurosci.</i> 2003; 20:173-88. Lindholm P, Voutilainen MH, Lauren J, et al. Novel neurotrophic factor CDNF protects and rescues midbrain dopamine neurons in vivo. <i>Nature</i> 2007; 448:73-7.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4349P; MANF Peptide Cat.No. PK-AB718-1463; Rat Brain Tissue Lysate

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