

# Neurotrypsin (NT) antibody (pAb)

## Rabbit Anti-Human/Mouse/Rat Neurotrypsin (NT)

### Instruction Manual

Catalog Number	PK-AB718-4105
Synonyms	Neurotrypsin Antibody: Motopsin, Protease serine 12, PRSS12, MRT1, BSSP3
Description	Neurotrypsin is a central nervous system-expressed serine protease whose truncation or absence causes nonsyndromic mental retardation. It is most prominently expressed in structures that are involved in the processing and storage of learned behavior and memory, such as the cerebral cortex, the hippocampus, and amygdala. Evidence suggests that neurotrypsin has multiple functions, including axonal outgrowth, maintaining neuronal plasticity, and arranging the perineuronal environment, partly in coordination with other proteases including tissue plasminogen activator. At least two isoforms of neurotrypsin are known to exist.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Neurotrypsin antibody was raised in rabbits against a 23 amino acid peptide from near the amino terminus of human neurotrypsin.
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.  Neurotrypsin antibody can be used for detection of neurotrypsin by Western blot at 0.5 - 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	Gschwend TP, Krueger SR, Kozlov SV, et al. Neurotrypsin, a novel multidomain serine protease expressed in the nervous system. <i>Mol. Cell. Neurosci.</i> 1997; 9:207-19. Molinari F, Rio M, Meskanaite V, et al. Truncating neurotrypsin mutation in autosomal recessive nonsyndromic mental retardation. <i>Science</i> 2002; 1779-81. Mitsui S, Yamaguchi N, Osako Y, et al. Enzymatic properties and localization of motopsin (PRSS12), a protease whose absence causes mental retardation. <i>Brain Res.</i> 2007; 1136:1-12.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4105P; Neurotrypsin Peptide Cat.No. PK-AB718-1220; SK-N-SH Cell Lysate

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