

# CAPS2 (IN) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat CAPS2 (Calcium-dependent secretion activator 2, CADPS2)

## Instruction Manual

Catalog Number	PK-AB718-4565
Synonyms	CAPS2 Antibody: Calcium-dependent secretion activator 2, CADPS2
Description	CAPS2 and the related protein CAPS1 encode novel neural/endocrine-specific cytosolic and peripheral membrane proteins. Both are essential components of the synaptic vesicle priming machinery and are required for the Ca <sup>2+</sup> -regulated exocytosis of secretory vesicles; CAPS-deficient neurons contain no or very few fusion competent synaptic vesicles, causing a selective impairment of fast phasic transmitter release. Like CAPS1, numerous isoforms of CAPS2 are known to exist. These isoforms show distinct expression patterns in the brain. For example, CAPS2b showed high expression in the developing cerebellum. Furthermore, one version of CAPS2 mRNA that lacks exon 3 is detected in some autistic patients, suggesting that the differential expression pattern of CAPS2 is involved in neuronal development. This CAPS2 antibody is predicted to be specific to CAPS2 and not recognize CAPS1.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	CAPS2 antibody was raised in rabbits against a 19 amino acid peptide near the center of the human CAPS2.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	This CAPS2 antibody is predicted to be specific to CAPS2 and not recognize CAPS1.
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 <b>INote:</b> Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually.  CAPS2 antibody can be used for detection of CAPS2 by Western blot at 0.5 - 1 µg/mL. Antibody can also be used for immunohistochemistry starting at 5 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Walent JH, Porter BW, and Martin TF. A novel 145 kD brain cytosolic protein reconstitutes Ca(2+)-regulated secretion in permeable neuroendocrine cells. Cell 1992; 70:765-775. Juckusch WJ, Speidel D, Sigler A, et al. CAPS-1 and CAPS-2 are essential synaptic vesicle priming proteins. Cell 2007; 131:796-808. Sadakata T, Washida M, and Furuichi T. Alternative splicing variants in mouse CAPS2: differential expression and functional properties of splicing variants. BMC Neurosci. 2007; 8:25-35.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4565P; CAPS2 Peptide Cat. No. PK-AB718-1303; Human Brain Tissue Lysate

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