

# SCRAPPER (NT) antibody (pAb)

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

### Instruction Manual

Catalog Number	PK-AB718-4451
Synonyms	SCRAPPER Antibody: F-box, Leucine rich repeat protein 20, FBXL20, FBI2, FBL20
Description	Members of the F-box protein family, such as Scrapper, are characterized by an approximately 40-amino acid F-box motif. SCF complexes, formed by SKP1, cullin, and F-box proteins, act as protein-ubiquitin ligases. Scrapper is selectively expressed in the brain, broadly expressed within the mouse CNS and is abundant at presynaptic membrane. Scrapper has orthologs in <i>C. elegans</i> , <i>D. melanogaster</i> , and mammals which suggests that it might function as an important membrane-localized E3 ligase in various species. Scrapper is a major presynaptic E3 ubiquitin ligase that acts through RIM1a via degradation and the ubiquitin-proteasome-system (UPS) pathway to critically regulate synaptic transmission. This identifies protein degradation as a mechanism for holding synaptic communication in check.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	SCRAPPER antibody was raised in rabbits against an 18 amino acid peptide near the amino terminus of the human SCRAPPER.
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 <b>INote:</b> Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. SCRAPPER antibody can be used for detection of SCRAPPER by Western blot at 0.5 - 1 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 MS, Tsai PI, and Chien CT. F-box proteins: the key to protein degradation. <i>J. Biomed. Sci.</i> 2006; 13:181-91. Dobie F and Craig AM. A fight for neurotransmission: SCRAPPER trashes RIM. <i>Cell</i> 2007; 130:775-7. Yao I, Takagi H, Ageta H, et al. SCRAPPER-dependent ubiquitination of active zone protein RIM1 regulates synaptic vesicle release. <i>Cell</i> 2007; 130:943-57. Ding M, Chao D, Wang G, et al. Spatial regulation of an E3 ubiquitin ligase directs selective synapse elimination. <i>Science</i> 2007; 317:947-51.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4451P; SCRAPPER Peptide Cat.No. PK-AB718-1288; A20 Cell Lysate

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