

SPAST monoclonal antibody

Other name: SPG4, ADPSP, FSP2

Cat. No. C15200001

Type: Monoclonal

Isotype: IgG2a

Source: Mouse

Lot #: 001

Size: 50 µg/50 µl

Concentration: 1 µg/µl

Specificity: Human: positive. Other species: not tested.

Purity: Affinity purified monoclonal antibody in PBS containing 0.05% azide.

Storage: Store at -20°C; for long storage, store at -80°C. Avoid multiple freeze-thaw cycles

Precautions: This product is for research use only. Not for use in diagnostic or therapeutic procedures

Description: Monoclonal antibody raised in mouse against SPAST (Spastin), using the full length recombinant protein.

Applications

	Suggested dilution	Results
Western blotting	1:500	Fig 1

Target description

SPAST (UniProt/Swiss-Prot entry Q9UBP0) is a member of the AAA (ATPases associated) protein family. Members of this protein family share an ATPase domain and have functions in a variety of cellular activities including membrane trafficking, intracellular motility, organelle biogenesis, protein folding, and proteolysis. SPAST seems to have a role in microtubule dynamics through its function as a microtubule-severing protein. Mutations in the ATPase binding domain of SPAST cause hereditary spastic paraplegias (HSP), a large group of clinically similar disorders.

Results

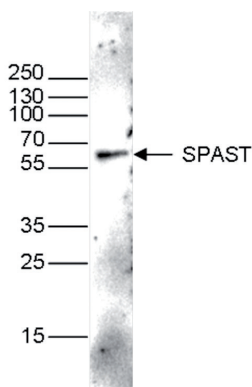


Figure 1. Western blot analysis using the Diagenode monoclonal antibody directed against SPAST

Whole cell extracts from HeLa cells were analysed by Western blot using the Diagenode antibody against SPAST (Cat. No. C15200001) diluted 1:500. The position of the protein of interest is indicated on the right; the marker (in kDa) is shown on the left.

Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK
Rue Bois Saint-Jean, 3
4102 Seraing (Ougrée) - Belgium
Tel: +32 4 364 20 50
Fax: +32 4 364 20 51
orders@diagenode.com
info@diagenode.com

Diagenode Inc. USA | NORTH AMERICA

400 Morris Avenue, Suite 101
Denville, NJ 07834 - USA
Tel: +1 862 209-4680
Fax: +1 862 209-4681
orders.na@diagenode.com
info.na@diagenode.com

Last update: September 3, 2015