

LC20 Protein

Full-length mouse recombinant protein expressed in E. coli cells

Catalog # M89-54G

Lot # V2598-4

Product Description

Recombinant full-length mouse LC20 was expressed in E.coli cells using an N-terminal GST tag. The gene accession number is [BC055439](#).

Gene Aliases

MYL9, MGC3505, MLC2, MRLC1, MYRL2

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

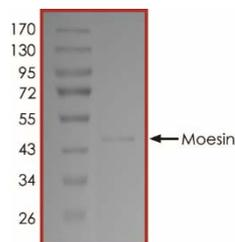
Scientific Background

LC20 is the myosin light chain that may regulate muscle contraction by modulating the ATPase activity of myosin heads (1). LC20 protein binds calcium and is activated by myosin light chain kinase. Two transcript variants encoding different isoforms have been found for LC20 and the deduced 172-amino acid protein is highly conserved, with only 3 differences between the human and chicken proteins. Light chain phosphorylation causes the folded monomeric form of myosin to extend and assemble into filaments. This observation established the involvement of the LC20 in conformational transitions of smooth muscle myosin (2).

References

1. Wagner, P D. et al: Regulation of the actin-activated ATPase of aorta smooth muscle myosin. J Biol Chem. 1986 Jun 15;261(17):7778-83.
2. Trybus, K M. et al: The regulatory light chain is required for folding of smooth muscle myosin. J Biol Chem. 1988 Nov 5;263(31):16485-92.

Purity



The purity was determined to be **>95%** by densitometry. Approx. MW **45kDa**.

LC20 Protein

Full-length mouse recombinant protein expressed in E. coli cells

Catalog #	M89-54G
Lot #	V2598-4
Purity	>95%
Concentration	0.5 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.