

MO25 α Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog # M51-30G

Lot # C125-1

Product Description

Recombinant full-length human MO25 α was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_016289](#).

Gene Aliases

CAB39, CGI-66, FLJ22682

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

MO25 α (mouse protein 25 alpha) is a 40-kDa protein that interacts with the STE20-related adaptor-alpha (STRAD α) pseudo kinase to form a regulatory complex capable of stimulating the activity of the LKB1 tumor suppressor protein kinase (1). LKB1 plays a critical role in cell proliferation, polarity and energy metabolism. LKB1 is mutated in the inherited Peutz-Jeghers cancer syndrome (PJS). MO25 α binds directly to a conserved Trp-Glu-Phe sequence at the STRAD α C terminus, and markedly enhances the binding of STRAD α to LKB1 thereby increasing LKB1 catalytic activity (2).

References

1. Boudeau, J. et al: Analysis of the LKB1-STRAD-MO25 complex. *J Cell Sci.* 2004;117(Pt 26):6365-75.
2. Milburn, C.C. et al: Crystal structure of MO25 alpha in complex with the C terminus of the pseudo kinase STE20-related adaptor. *Nat. Struct. Mol. Biol.* 2004;11(2):193-200.

Purity



The purity of MO25 α was determined to be **>95%** by densitometry. Approx. MW **63kDa**.

MO25 α Protein

Full-length recombinant protein expressed in Sf9 cells

Catalog Number M51-30G

Specific Lot Number C125-1

| | |
|--------------------|--|
| Purity | >95% |
| Concentration | 0.2 µg/µl |
| Stability | 1 yr 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.