

MBP Protein

Full length recombinant human protein expressed E. coli cells

Catalog # M42-54G

Lot # Q2560-7

Product Description

Recombinant full length human MBP was expressed in E.coli cells using an N-terminal GST tag. The gene accession number is [NM_001025090](#).

Gene Aliases

None

Formulation

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

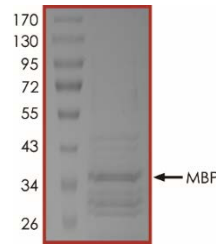
Scientific Background

MBP or Myelin Basic Protein forms dimers across the cytoplasmic apposition during the formation of myelin. The region of the protein involved in folding, polymerization and substrate specificities is conserved in various species and it may have a specialized role in protein-lipid interactions in the myelin membrane (1). MBP is an efficient substrate for numerous protein kinases and share similarities with the MARCKS protein in terms of having extended conformations regulated by their environment, N-terminal modifications, a dual nature of interactions with lipids, and binding to actin and Ca²⁺-calmodulin (2).

References

1. Chapman, B.E. et al: Conformation of myelin basic protein and its role in myelin formation. Adv Exp Med Biol. 1978;100:207-20.
2. Harauz, G. et al: Analogous structural motifs in myelin basic protein and in MARCKS. Mol Cell Biochem.2000 Jun;209(1-2):155-63.

Purity



The purity was determined to be **>65%** by densitometry. Approx. MW **36kDa**.

MBP Protein

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

| | |
|--------------------|---|
| Catalog # | M42-54G |
| Lot # | Q2560-7 |
| Purity | >65% |
| Concentration | 0.1 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.