

Myelin basic protein (human brain), (high purity)

MBP is a peripheral membrane protein, and its interaction with lipids is generally believed to be critical for the formation and stability of the multilamellar myelin sheath. It has been identified as a Ca^{2+} -calmodulin regulated agent for actin polymerization, tropomyosin and actomyosin function, tubulin stabilization, and clathrin assembly. MBP is generally considered to be the antigen responsible for autoimmunity in multiple sclerosis (MS) and it has the capability to induce experimental allergic encephalomyelitis (EAE). MBP is also an endogenous inhibitor of the high-affinity cannabinoid binding site in the brain and a substrate for numerous protein kinases (MAPK, serine/threonine kinase, protein kinase C). MBP is also suggested to be involved in intracellular signalling and as cytoskeletal components.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-200-606-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for 12 months at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	MBP
Appearance	White to off-white powder.
Application Notes	Can be used to induce relapsing experimental autoimmune encephalomyelitis in animals. This product may also be used as a substrate for protein kinases. Suitable for use in ELISA or for antigen applications in immunological protocols. Immunoreactivity is verified using polyclonal and monoclonal antibodies against MBP.
Concentration	1mg/ml after reconstitution.
Formulation	Lyophilized. No preservatives added.
MW	MBP contains 14.2kDa (13.0 minor) and 18.5kDa (17.2 minor) isoforms as well as their dimers (13.5% SDS-PAGE, non-reducing conditions).
Purity	≥95% (SDS-PAGE; Immunoblot)
Reconstitution	Reconstitute with 10mM HCl. Further dilutions can be made in assay buffer.
Source	Purified from human brain.
UniProt ID	P02686



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)