

## Data Sheet

### MBP

Human, recombinant, N-terminal His-tag

**Catalog #:** 40535

**Lot #:** 130517-Q2      **Conc.:** 1.1 mg/ml

**Formulated in:** 40 mM HEPES pH 7.5, 250 mM NaCl, and 20% glycerol.

**Stability:** >6 months at -70°C. *Avoid multiple freeze/thaw cycles.*

#### References:

1. Kishimoto, A., *et al.*, *J. Biol. Chem.* 1985;**260(23)**:12492-12499
2. Quandt, J.A., *et al.*, *J. Immunol.* 2012;**189(6)**:2897-2908

#### Description:

Human myelin basic protein (MBP), GenBank Accession No. NM\_001025101, a.a. 2-304(end), with N-terminal His-tag, MW=34 kDa, expressed in an *E. coli* cell expression system.

#### Application:

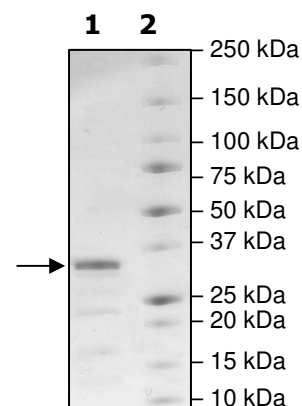
Useful as an immunogen or as a substrate for phosphorylation, citrullination, or methylation studies.

### Quality Assurance

#### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
2 µg MBP  
**Lane 2:**  
Protein Marker

**MW:** 34 kDa  
**Purity:** ≥70%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**  
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6044 Cornerstone Court West, Suite E  
San Diego, CA 92121  
Tel: 1.858.829.3082  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**Sequence:**

**MHHHHHHHGNHAGKRELNAEKASTNSETNRGESEKKRNLGELSRTTSEDNEVFGEADA  
NQNGTSSQDTAVTDSKRTADPKNAWQDAHPADPGSRPHLIRLFSRDAPGREDNTFK  
DRPSEDELQTIQEDSAATSESLDVMASQKRPSQRHGSKYLATASTMDHARHGFLPRH  
RDTGILDSIGRFFGGDRGAPKRGSGKDSHHPARTAHYGSLPQKSHGRTQDENPVVHFF  
KNIVTPRTPPPSQGKGRGLSLSRFSWGAEGQRPGFGYGGGRASDYKSAHKGFKGVDAQ  
GTLKIFKLGGRDSRSGSPMARR**

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**  
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)