

Data Sheet

EIF4EBP1

Human, recombinant, N-terminal His-tag

Catalog #: 40510

Lot#: 120103 **Conc.:** 0.22 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCL, 200 mM
imidazole, 20% Glycerol and 3mM DTT.

Stability: >12 months at -80°C

References:

1. Fonseca, B.D. et al., *J. Biol. Chem.* **282** (34), 24514-24524 (2007).
2. Armengol, G. et al., *Cancer Res.* **67** (16), 7551-7555 (2007).

Description:

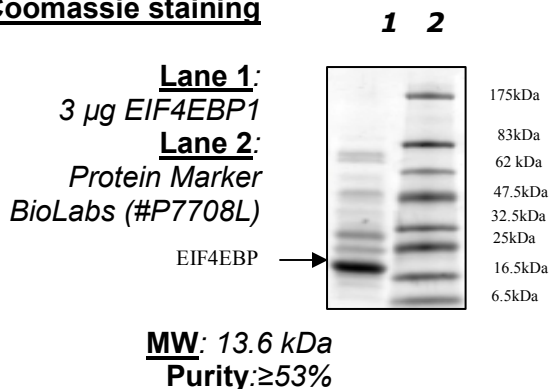
Human EIF4EBP1 (eukaryotic translation initiation factor 4E binding protein 1) (GenBank Accession No. NM_004095), full length with N-terminal His-tag, MW= 13.6 kDa, expressed in an E. Coli expression system.

Application:

Suitable substrate for mTOR

Quality Assurance

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



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 1.858.259.5734

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com