

Data Sheet

Elob, FLAG-Tag

Human, Recombinant, N-terminal FLAG-tag

Catalog #: 100256

Lot #: 190321

Conc.: 0.24 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 80 µg/ml FLAG peptide, 3mM DTT.

Description: Human Elob also known as Elongin B, RNA Polymerase II Transcription Factor SIII Subunit B, TCEB2, and SIII, GenBank Accession No. NM_007108, a.a. 2-118(end), with N-terminal FLAG-tag, expressed in a HEK293 expression system, MW=14 kDa.

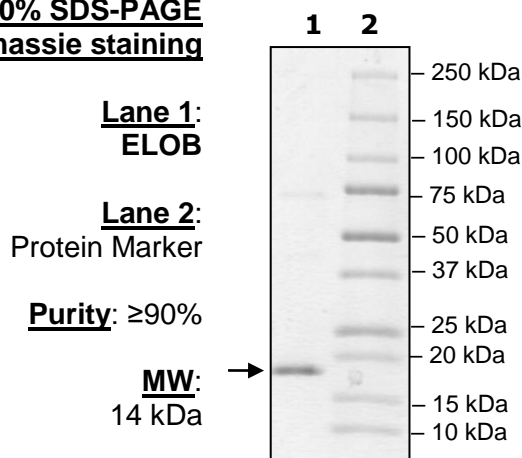
Stability: At least 6 months at -80°C.
Avoid freeze/thaw cycles.

References:

1. Kibel, A. *et al.*, *Science*. **269(5229)** : 1444-1446 (1995).
2. Ohh, M. *et al.*, *JCI*. **104**: 1583-1591 (1999).

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 **1.858.202.1401** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com