

Data Sheet

SOS1, FLAG-Tag, Avi-Tag

Human, Recombinant, N-terminal FLAG-tag,
C-terminal Avi-Tag

Catalog #: 100752

Lot #: 200828

Conc.: 0.55 mg/ml

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

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

References:

1. Shi, H., *et al.*, *Plant Cell*. 2002; **14**:
465-477.
2. Shi, H., *et al.*, *PNAS*. 2000; **97(12)**:
6896-6901.

Description: Human Son of sevenless
homolog 1 (SOS1), also known as guanine
nucleotide exchange factor 1 (GF1),
GenBank Accession No. NM_005633, a.a.
594-1049, with N-terminal FLAG-tag and
C-terminal Avi-tag, expressed in a
Baculovirus infected Sf9 cell expression
system, MW=57 kDa.

Quality Assurance

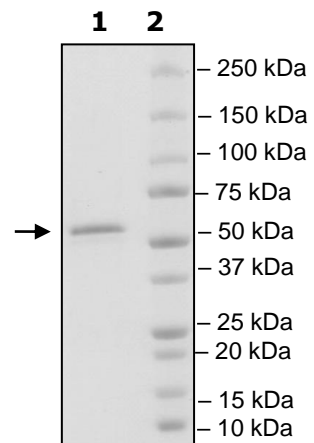
4-20% SDS-PAGE **Coomassie staining**

Lane 1:
SOS1

Lane 2:
Protein Marker →

Purity: >90%

MW: 57 kDa



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 **1.858.259.5734**

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

Please visit our website at: www.bpsbioscience.com