

## Data Sheet

### SOS1, FLAG-Tag, Avi-Tag, Biotin-Labeled

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

**Catalog #:** 100753

**Lot #:** 200602

**Conc.:** 1.31 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. This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%.

## Quality Assurance

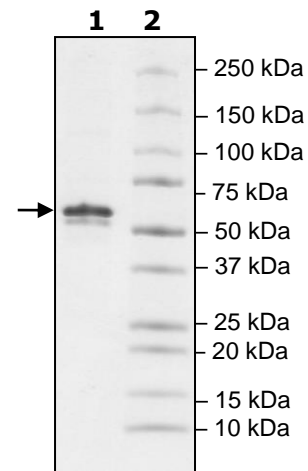
### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
**SOS1**

**Lane 2:**  
Protein Marker

**Purity:** >85%

**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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)