

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

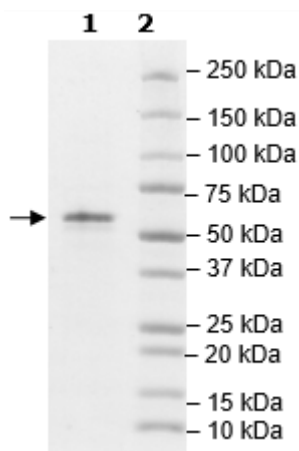
Catalog: 100753  
Lot: 200603

## Product Information

<b>Description:</b>	Human Son of sevenless homolog 1 (SOS1), also known as guanine nucleotide exchange factor 1 (GEF1), encompassing amino acids 594-1049. This construct contains an N-terminal FLAG-tag and a C-terminal Avi-tag™. The recombinant protein was affinity purified and enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%.
<b>Species:</b>	Human
<b>Construct:</b>	SOS1 (FLAG-594-1049-Avi)-(Biotin)
<b>Concentration:</b>	0.59 mg/ml
<b>Expression System:</b>	Sf9
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 3 mM DTT, and 80 µg/ml FLAG peptide
<b>MW:</b>	57 kDa
<b>Genbank Accession:</b>	NM_005633
<b>Label:</b>	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation confirmed to be ≥90%.
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.

## Quality Control Data

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



### Biotin-Avidin Pulldown

