

Data Sheet

SARM1, FLAG-Tag

Human, Recombinant, N-terminal FLAG-Tag

Catalog #: 100069

Lot #: 180723 **Conc.:** 0.56 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 3mM DTT, 100ug/ml FLAG peptide

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

References:

1. Osterloh, J. *et al.*, *Science*. Jul; **337(6093)**: 481-484 (2012).
2. Molday, L. *et al.*, *JBC*. **282**: 32792-32801. (2007).

Description: Human SARM1, also known as Sterile Alpha and TIR Motif Containing 1, and SAMD2, a.a. 28-724(end), with N-terminal FLAG-tag, GenBank Accession No. NM_015077, MW = 77 kDa, expressed in a HEK293 expression system.

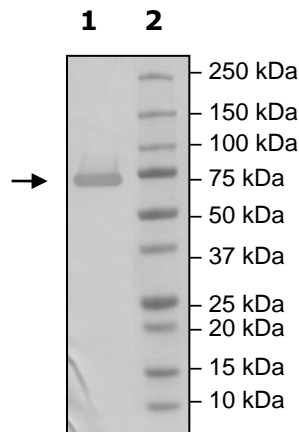
Application: Useful as a positive control for Western blot studies.

Quality Assurance

Western Blot
w/ FLAG-anti body

Lane 1:
SARM1
Lane 2:
Protein Marker

MW: 77 kDa

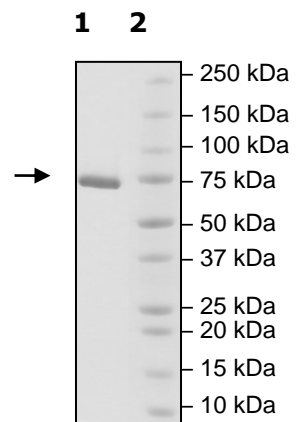


4-20% SDS-PAGE
Coomassie staining

Lane 1:
SARM1
Lane 2:
Protein Marker

Purity: ≥90%

MW: 77 kDa



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com