

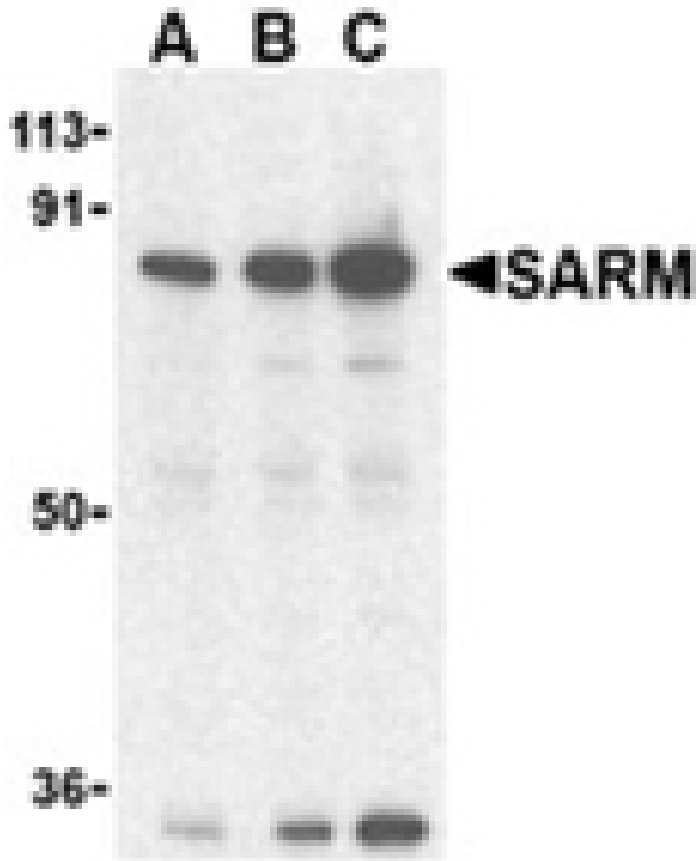
SARM polyclonal antibody

SARM (SAM and ARM-containing protein) is thought to serve as an adaptor protein for the Toll-like receptors (TLRs) and allow for the activation of downstream kinases and NF- κ B, ultimately leading to the expression of proteins involved in host defense. While SARM has not been conclusively shown to associate directly with TLRs, the presence of a Toll-interleukin-1 (TIR) domain in SARM is consistent with a role as a signaling molecule.

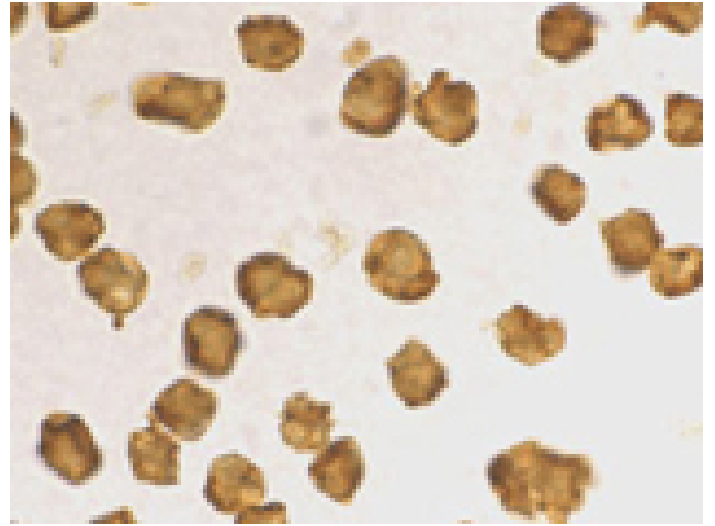
This antibody is covered by our [Worry-Free Guarantee](#).

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of SARM in Daudi cell lysates with SARM antibody at (A) 0.5, (B) 1, and (C) 2 μ g/ml.



Immunocytochemistry analysis of Daudi cells using SARM antibody at 2 μ g/ml.

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	TIR-1, Sterile α and TIR motif-containing protein 1, SAM and ARM-containing protein, Sterile α motif and armadillo repeat-containing protein
Application	ICC, WB
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Gene/Protein Identifier	NP_055892 (RefSeq)
Host	Rabbit
Immunogen	Synthetic peptide corresponding to the sequence near the C-terminus of human SARM.
Purity Detail	Ion exchange purified.
Recommendation Dilutions/Conditions	Immunocytochemistry (2 μ g/ml)Western Blot (0.5-2.0 μ g/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse
UniProt ID	Q6SZW1
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024
ENZO LIFE SCIENCES
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com