

SARM (human) monoclonal antibody (Sarmy-1)

In mammals, TIR domain adapters, including Myd88, TIRAP, TICAM-1 (also called TRIF), and TRAM, act in Toll-like receptor (TLR) signaling pathway. The last member of this family called SARM (sterile and armadillo motifs) is highly conserved but has no known function in any organism. SARM is the only mammalian TIR-domain adapter proteins conserved in *C. elegans*, suggesting that SARM is the most ancient member of this class of signaling proteins. Overexpression of SARM fails to activate either NF- κ B- or IRF-3-dependent reporter gene expression, two transcription factors that are sensitive to TLR signaling, suggesting that SARM is functionally distinct from other mammalian TIR domain-containing proteins.

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

Ordering Information

[Order Online »](#)

ALX-804-831-C100	100 μ g
------------------	-------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored at -20°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°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	IP
Application Notes	Does not work for Western Blot.
Clone	Sarmy-1
Formulation	Liquid. In PBS containing 10% glycerol and 0.02% sodium azide.
Host	Mouse
Immunogen	Recombinant human SARM (aa 93-292).
Isotype	IgG2b
Recommendation Dilutions/Conditions	Immunoprecipitation (1:200)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from concentrated hybridoma tissue culture supernatant.
Species Reactivity	Human



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
Q6SZW1
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee.

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.