

TLR9 polyclonal antibody

The TLRs (Toll-like receptors) are a group of evolutionarily conserved pattern recognition molecules that play a key role in host defense during microbial infection by regulating both innate and adaptive immune responses. TLRs are type I transmembrane proteins with ectodomains containing interspersed leucine-rich repeat (LRR) motifs that are involved in the recognition of various microbial ligands.

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

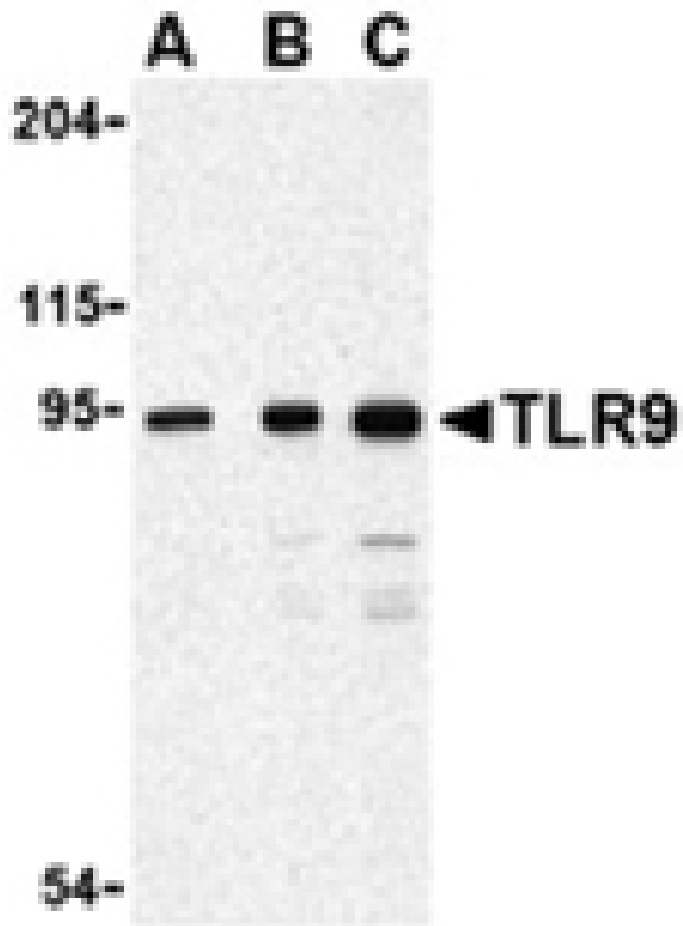
Ordering Information

[Order Online »](#)

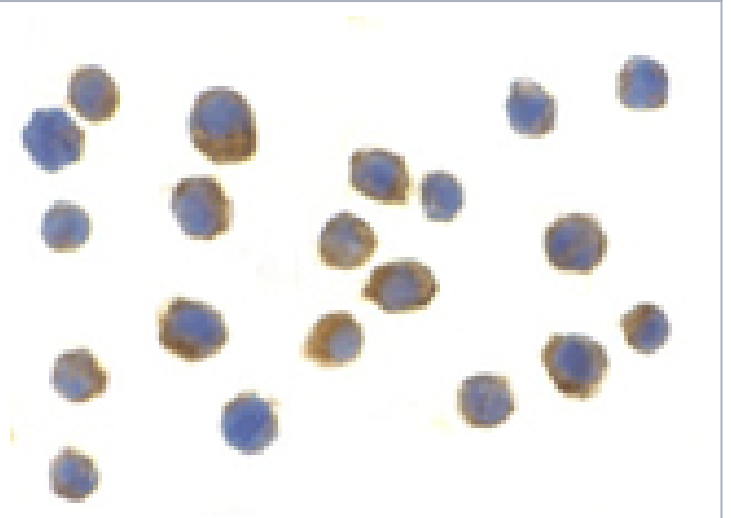
ADI-905-730-100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of TLR9 in Jurkat cell lysate with TLR9 antibody at (A) 0.5, (B) 1 and (C) 2µg/ml.



Immunocytochemistry analysis of TLR9 in Jurkat cells with TLR9 antibody at 2µg/ml.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Toll-like receptor 9
Application	ICC, WB
Application Notes	Detects a band of ~94kDa by Western blot.
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to the sequence near the middle of human TLR9.
Purity Detail	Peptide affinity purified.
Recommendation Dilutions/Conditions	Immunocytochemistry (2µg/ml)Western Blot (0.5-1.0µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse, Rat
UniProt ID	Q9NR96
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](mailto:info-usa@enzolifesciences.com)

European Sales Office
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-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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

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