

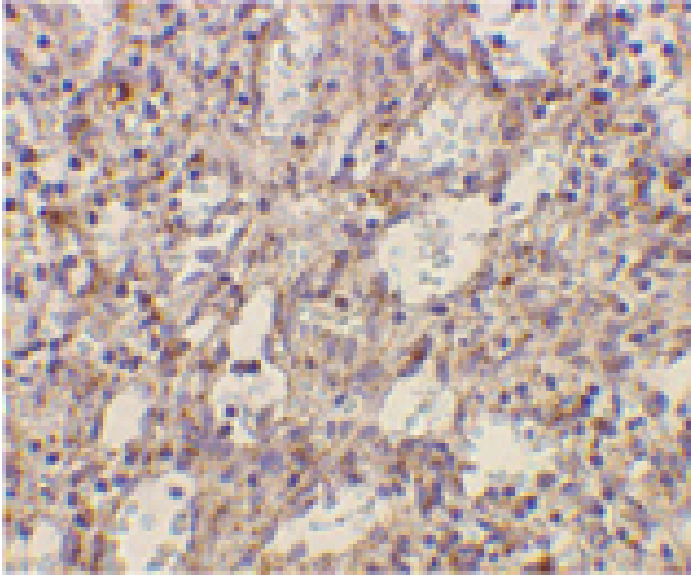
TLR2 polyclonal antibody

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.

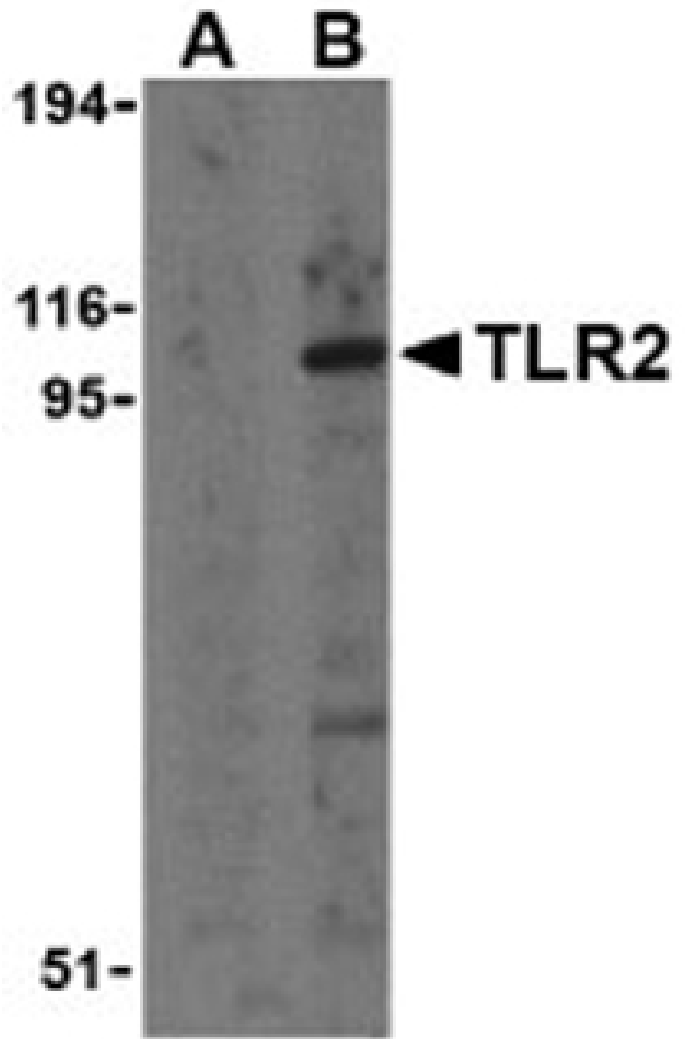
This antibody is covered by our [Worry-Free Guarantee](#).

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis of human spleen cells using TLR2 antibody at 2 μ g/ml.



Western blot analysis of TLR2 in A20 cell lysates with TLR2 antibody at 1 μ g/ml in the presence (A) and absence (B) of its blocking peptide.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Toll-like receptor 2
Application	IHC, WB
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to the sequence near the N-terminus of human TLR2.
Purity Detail	Ion exchange purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (2µg/ml)Western Blot (0.5-2.0µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse
UniProt ID	O60603
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-fr@enzolifesciences.com

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

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

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