

TLR7 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](#).

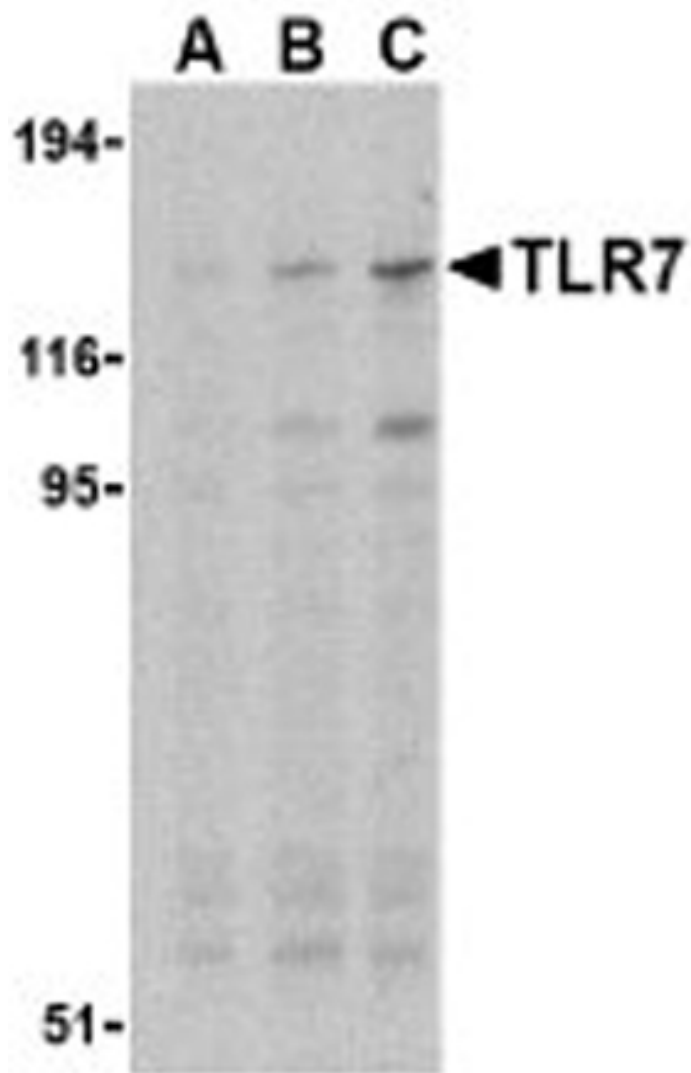
Ordering Information

[Order Online »](#)

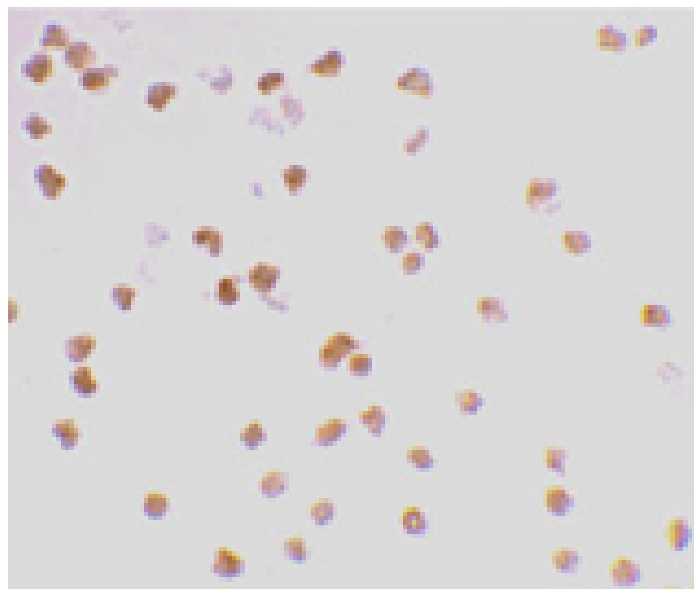
ADI-CSA-824-E	100µg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of TLR7 in Daudi cell lysates with TLR7 antibody at (A) 0.5, (B) 1, and (C) 2 μ g/ml.



Immunocytochemistry analysis of Daudi cells using TLR7 antibody at 2 μ g/ml.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Toll-like receptor 7

Application ICC, WB

Application Notes Detects a band of ~121kDa by Western blot.

Formulation Liquid. In PBS containing 0.02% sodium azide.

Gene/Protein Identifier NP_573474 (RefSeq)

Host Rabbit

Immunogen Synthetic peptide corresponding to the sequence near the middle of mouse TLR7.

Purity Detail Ion exchange purified.

Recommendation Immunocytochemistry (2µg/ml)Western Blot (0.5-2.0µg/ml)Suggested
Dilutions/Conditions dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human, Mouse

UniProt ID P58681

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



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-
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)