

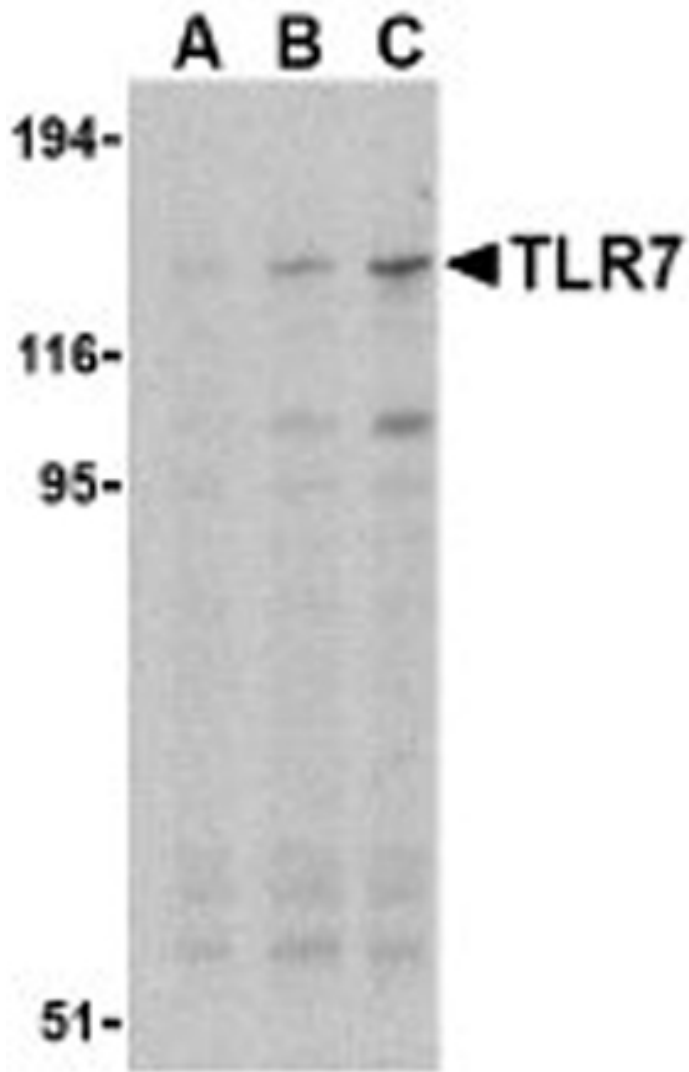
# TLR7 polyclonal antibody

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.

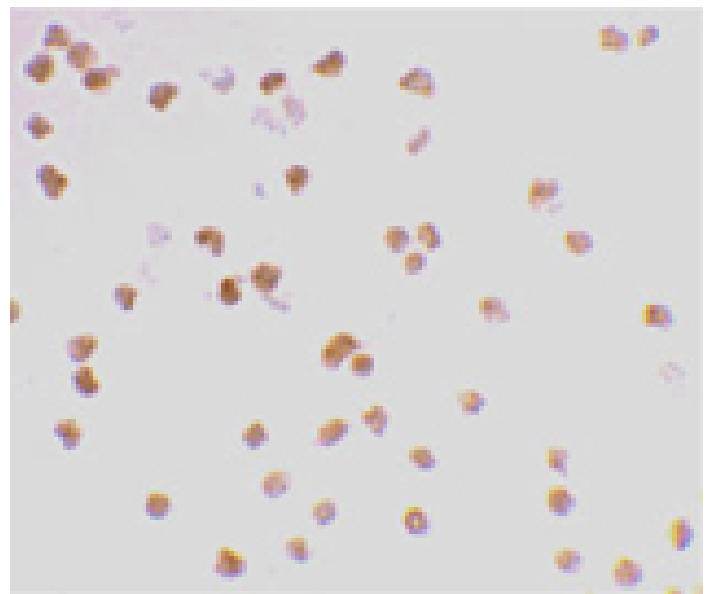
This antibody is covered by our [Worry-Free Guarantee](#).

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 $\mu$ g/ml.



Immunocytochemistry analysis of Daudi cells using TLR7 antibody at 2 $\mu$ 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 Dilutions/Conditions	Immunocytochemistry (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	P58681
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>