

Toll-like receptor 4 polyclonal antibody

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

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-210-638-C200	200µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

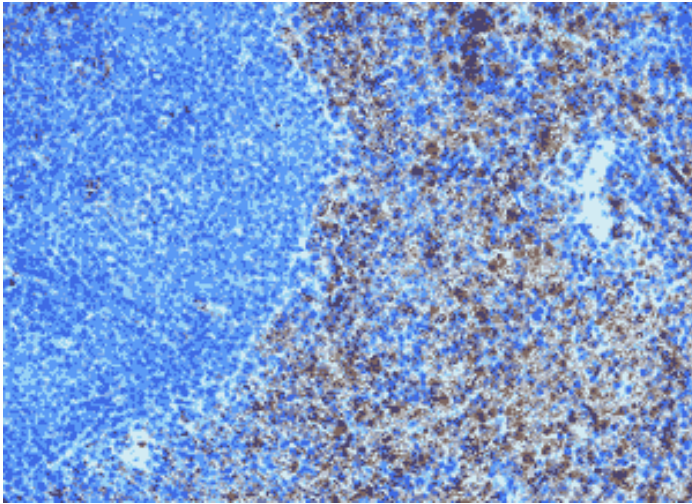


Figure 2: Immunohistochemistry of mouse TLR4 on paraffin section of mouse spleen using to Toll-like Receptor 4 (Prod. No. [ALX-210-638](#)).""",Toll-like receptor 4 polyclonal antibody
 Immunohistochemistry,Toll-like receptor 4 polyclonal antibody Immunohistochemistry,yes
 ALX-210-641,210-641.gif,<http://static.enzolifesciences.com/fileadmin/files/image/210-641.gif>,**Figure:** Immunohistochemistry on paraffin sections of mouse spleen using PAb to CCR3 (mouse) (Prod. No. [ALX-210-641](#)).""""""

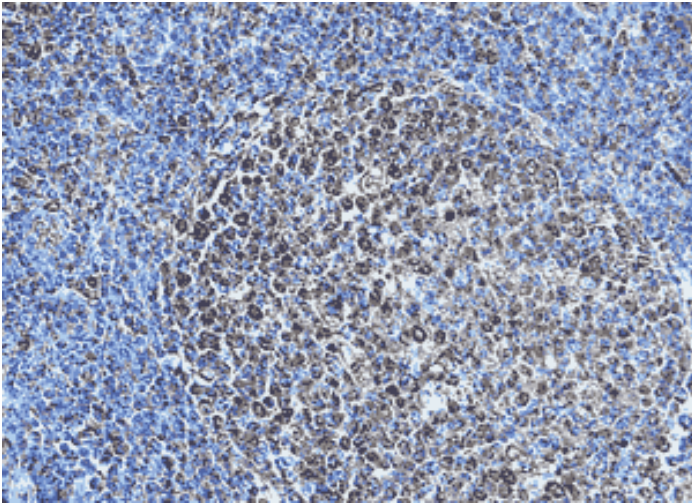


Figure 1: Immunohistochemistry of human TLR4 on paraffin section of human tonsil using to Toll-like Receptor 4 (Prod. No. ALX-210-638).

Handling & Storage

Handling	For maximum product recovery after thawing, centrifuge the vial before opening the cap. Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	CD284, TLR4
Application	ICC, IHC (PS), WB
Formulation	Liquid. In PBS containing 1mg/ml BSA and 0.1% sodium azide.
Host	Goat
Immunogen	Synthetic peptide corresponding to aa 161-192 (L161IQSFKLPEYFSNLTNLEHLDLSSNKIQSIYC192) of the N-terminal domain of human TLR4 (Toll-like receptor 4).
Purity Detail	Epitope-affinity purified IgG.
Recommendation Dilutions/Conditions	Immunocytochemistry (acetone fixed cells, >1:400)Immunohistochemistry (paraffin sections, >1:250)Western Blot (>1:1,000)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	O00206
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)