

Toll-like receptor 7 polyclonal antibody

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

Citations: 11

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

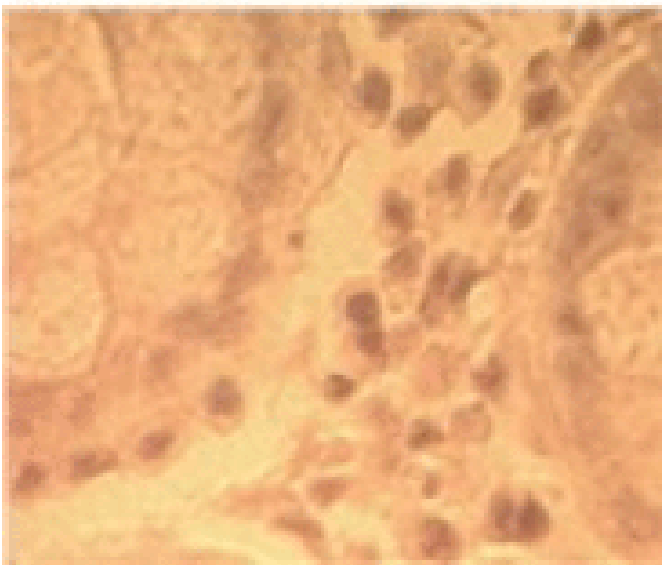
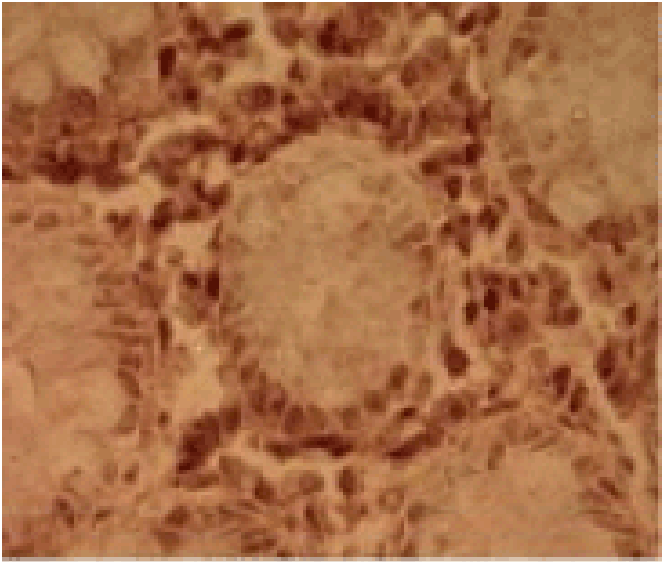


Figure 3: Immunohistochemical analysis of human colon tissue using PAb to Toll-like Receptor 7 (Prod. No. [ALX-210-874](#)) and an isotype control (bottom) at 5µg/ml. Toll-like receptor 7 polyclonal antibody Immunohistochemistry, Toll-like receptor 7 polyclonal antibody Immunohistochemistry, yes ALX-210-874, 210-

874_1.gif, http://static.enzolifesciences.com/fileadmin/files/image/210-874_1.gif, **Figure 1:** Western blot analysis of TLR7 in Raw cell lysate using Prod. No. [ALX-210-874](#) at 1µg/ml.

“””””””

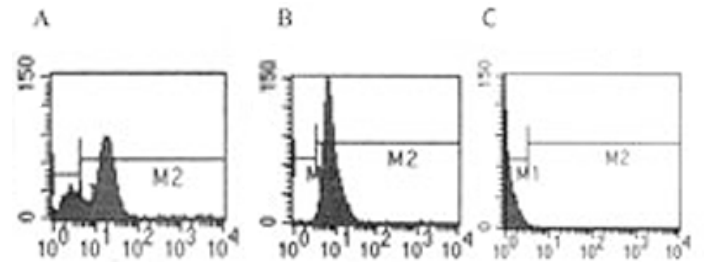


Figure 2: Intracellular staining by FACS analysis of TLR7 in Ramos cells (A) and Raw cells (B) using Prod. No. [ALX-210-874](#) at 2µg/ml. Histogram C shows the rabbit IgG isotype control. Toll-like receptor 7 polyclonal antibody Flow Cytometry, Toll-like receptor 7 polyclonal antibody Flow Cytometry, yes ALX-210-898, 210-

898.gif, <http://static.enzolifesciences.com/fileadmin/files/image/210-898.gif>, **Figure:** Western blot comparing three different PABs to VASP using 25µg human platelet protein. **Lane 1:** PAb to VASP (human) (Prod. N. ALX-210-898). **Lane 2:** PAb to VASP (M4) (Prod. N. ALX-210-725) (**Not available anymore!**). **Lane 3:** PAb to VASP (affinity purified) (Prod. N. ALX-210-880)., VASP (human) polyclonal antibody Western blot, VASP (human) polyclonal antibody Western blot, yes ALX-210-904, 210-

904.gif, <http://static.enzolifesciences.com/fileadmin/files/image/210-904.gif>, **Figure:** Western blot analysis of human and mouse cell lines with the anti-human NLRP1/NALP1 antibody (AL176) (Prod. No. [ALX-210-904](#)). **Method:** Total protein extracts (20µg) from various human (293-T

904.gif, <http://static.enzolifesciences.com/fileadmin/files/image/210-904.gif>, **Figure:** Western blot analysis of human and mouse cell lines with the anti-human NLRP1/NALP1 antibody (AL176) (Prod. No. [ALX-210-904](#)). **Method:** Total protein extracts (20µg) from various human (293-T

904.gif, <http://static.enzolifesciences.com/fileadmin/files/image/210-904.gif>, **Figure:** Western blot analysis of human and mouse cell lines with the anti-human NLRP1/NALP1 antibody (AL176) (Prod. No. [ALX-210-904](#)). **Method:** Total protein extracts (20µg) from various human (293-T

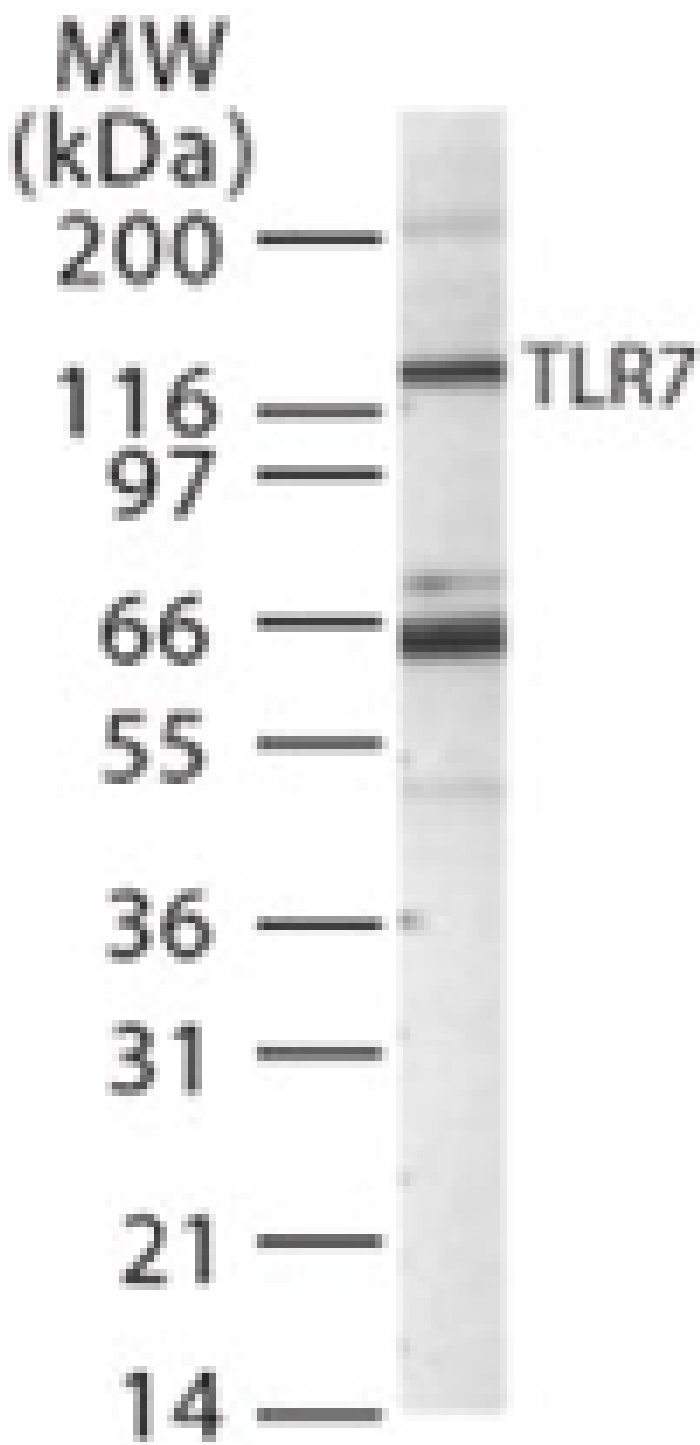


Figure 1: Western blot analysis of TLR7 in Raw cell lysate using Prod. No. ALX-210-874 at 1 μ g/ml.

Handling & Storage

Use/Stability	Stable for 6 months when stored at +4°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	TLR7
Application	Flow Cytometry, ICC, IHC (FS), IHC (PS), IP, WB
Application Notes	Flow Cytometry: suitable for cell surface and intracellular FC. Western Blot: in human Ramos cells, additional bands are seen
Formulation	Liquid. In PBS containing 0.05% BSA and 0.05% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 706-728 of human TLR7 (Toll-like receptor 7).
Purity Detail	Protein G-affinity purified.
Recommendation Dilutions/Conditions	Flow Cytometry (intracellular, 0.5-5µg/106 cells)Immunohistochemistry (paraffin sections, 5-10µg/ml)Western Blot (1-3µg/ml)
Species Reactivity	Human, Mouse, Rat
UniProt ID	Q9NYK1

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-
be@enzolifesciences.com](mailto:info-be@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)