

Toll-like receptor 9 monoclonal antibody (26C593.2)

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

Citations: 13

[View Online »](#)

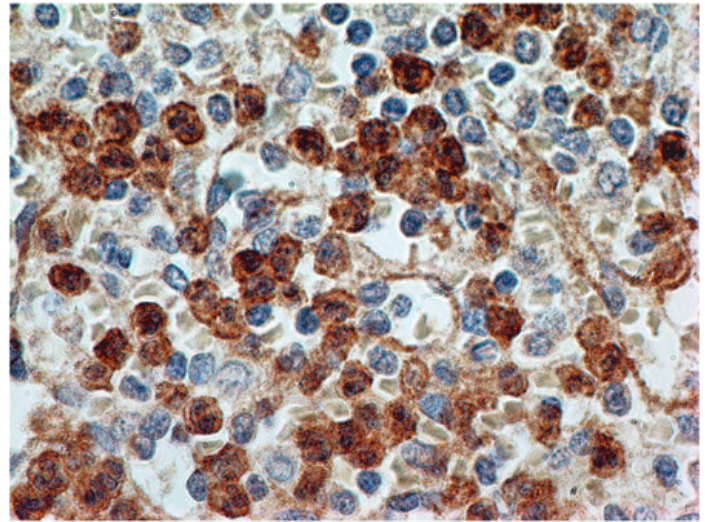
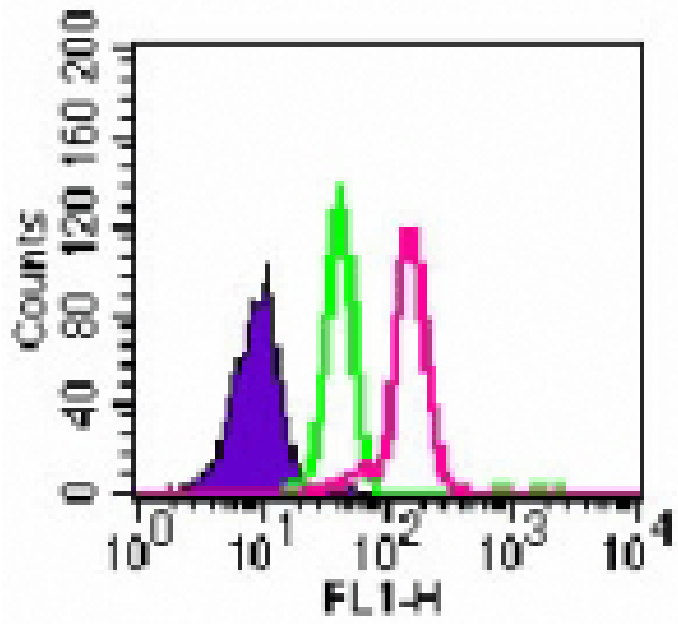
Ordering Information

[Order Online »](#)

ALX-804-364-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis: Formalin-fixed, paraffin-embedded human spleen probed with Toll-like receptor 9, mAb (26C593.2) at $5\mu\text{g/ml}$.

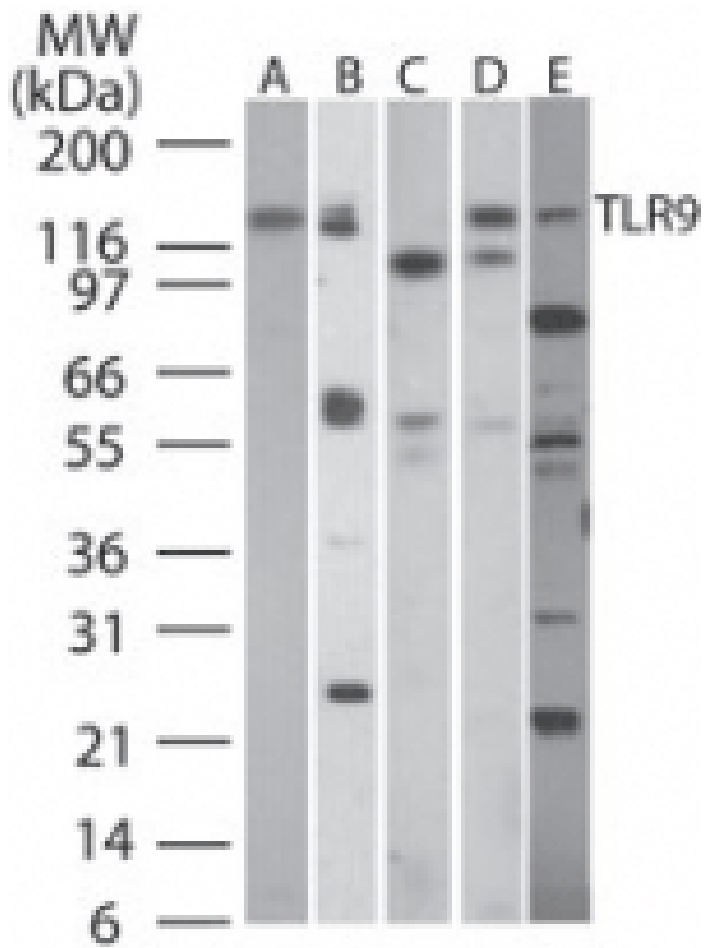


Figure 1: Western blot analysis of TLR9 using MAb to Toll-like Receptor 9 (human) (26C593.2) (Prod. No. ALX-804-364) at a dilution of 3 μ g/ml. A) human PBMC; B) human intestine; C) mouse intestine; D) rat intestine lysate; E) human MCF7 cell lysate (MAb dilution 5 μ g/ml).

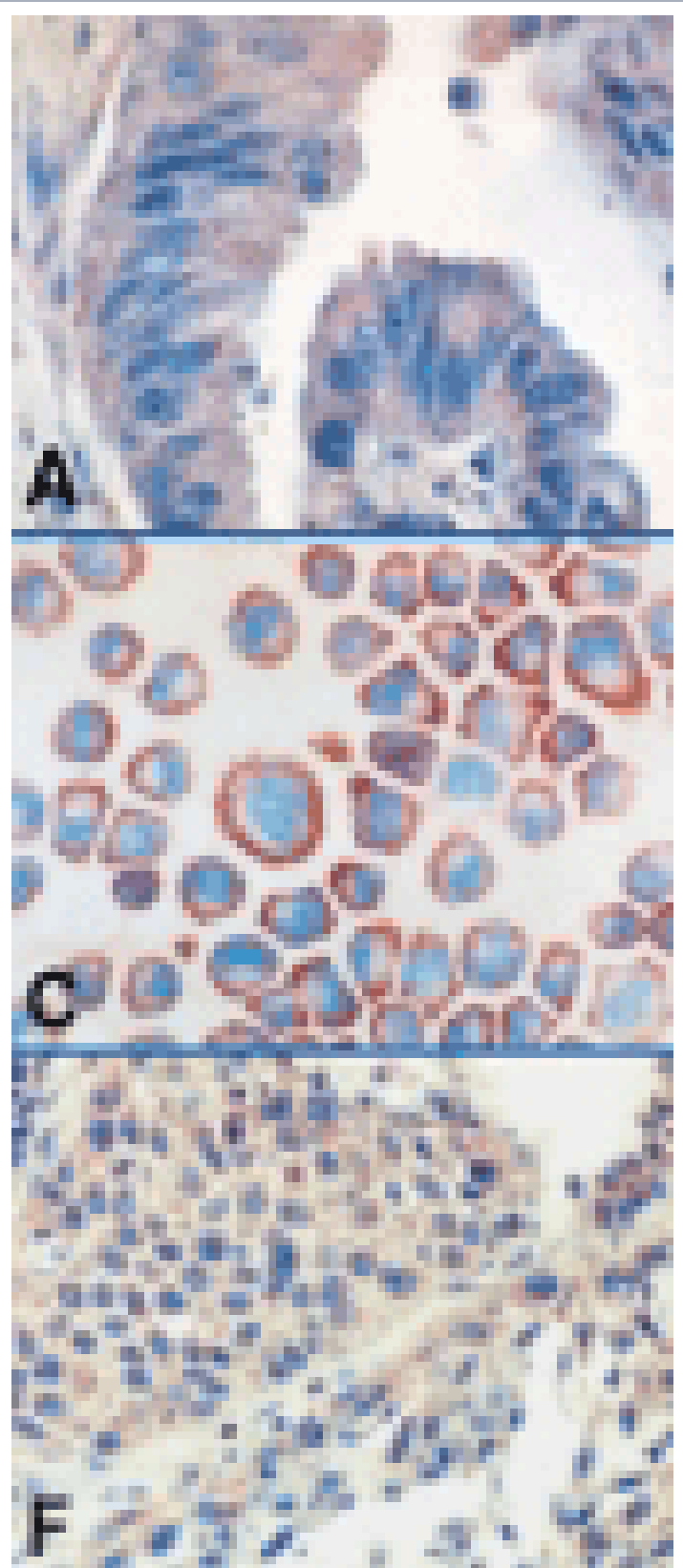


Figure 3: Immunohistochemical analysis of TLR9 in A) adenocarcinoma of the lung, C) A549 cells, and F) malignant lung tissues using MAb to Toll-like Receptor 9 (human) (26C593.2) (Prod. No. ALX-804-364) at 1:100 dilution, (600x magnification).

Handling & Storage

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	TLR9, CD289
Application	Flow Cytometry, IHC (FS), IHC (PS), WB
Application Notes	Flow Cytometry: Suitable for cell surface and intracellular FC. Detects a band of ~120kDa by Western blot, may also detect a smaller band of TLR9 isoform B.
Clone	26C593.2
Formulation	Liquid. In PBS containing 0.05% BSA and 0.05% sodium azide.
Host	Mouse
Immunogen	Synthetic peptide corresponding to aa 268-284 in the cytoplasmic portion of human TLR9 (Toll-like Receptor 9).
Isotype	IgG1
Positive Control	Whole cell lysate from human peripheral blood monocyte cells (PBMC) or Daudi can be used as positive control.
Purity Detail	Protein G-affinity purified.

Recommendation Dilutions/Conditions

Flow Cytometry (intracellular, 0.1-2µg/106 cells)Immunohistochemistry (frozen sections, 10-20µg/ml)Western Blot (1-5µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity

Dog, Horse, Human, Monkey, Mouse, Rat

UniProt ID

Q9NR96

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](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)