

Zap-70 monoclonal antibody (11H40)

Zeta-associated protein 70 (Zap-70) is a cytoplasmic protein tyrosine required for T cell antigen receptor (TCR) signaling and development. ZAP-70 interacts with the TCR by binding to tyrosine-phosphorylated immunoreceptor tyrosine-based activation motifs (ITAMs) present in the invariant subunits of the TCR complex.

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

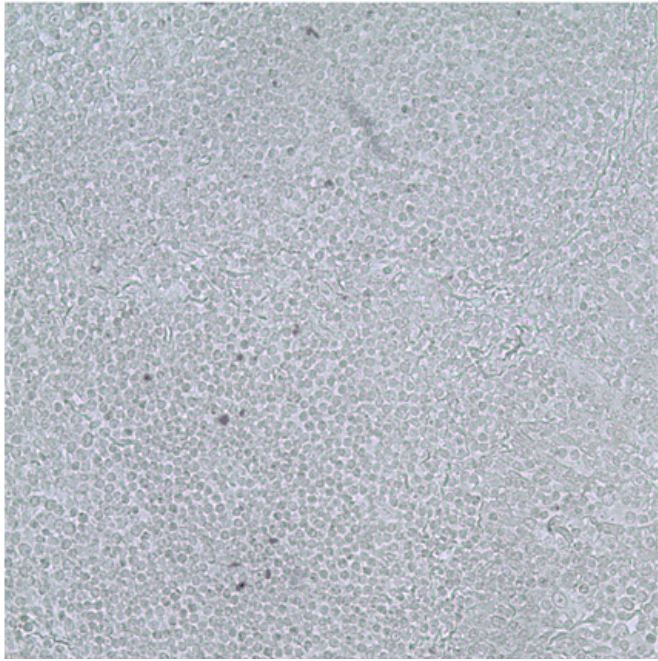
Ordering Information

[Order Online »](#)

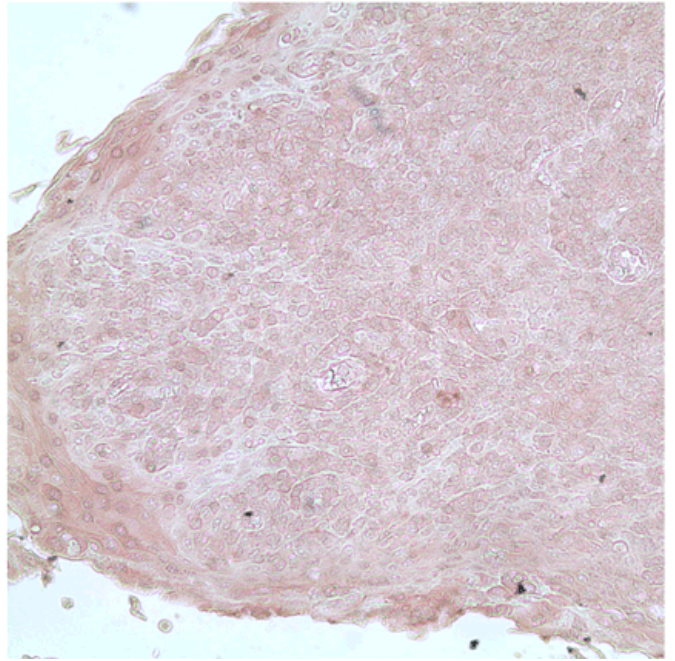
ENZ-ABS709-0100	100µg
-----------------	-------

Manuals, SDS & CofA

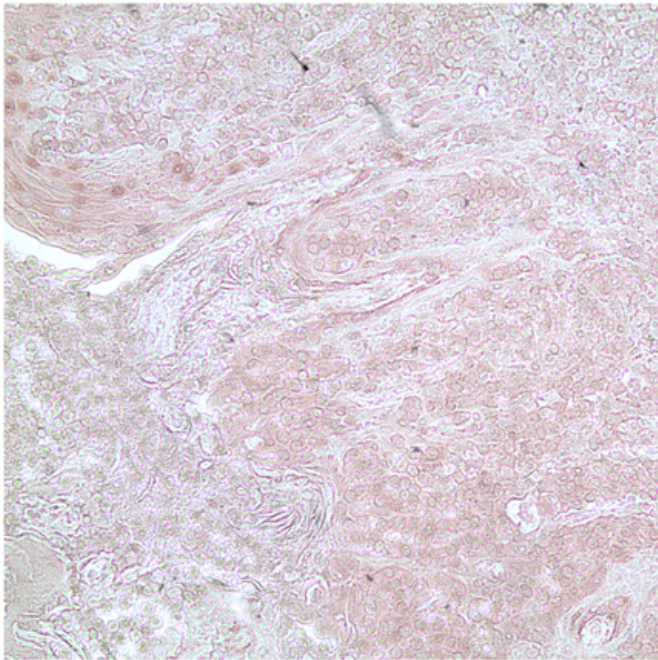
[View Online »](#)

A

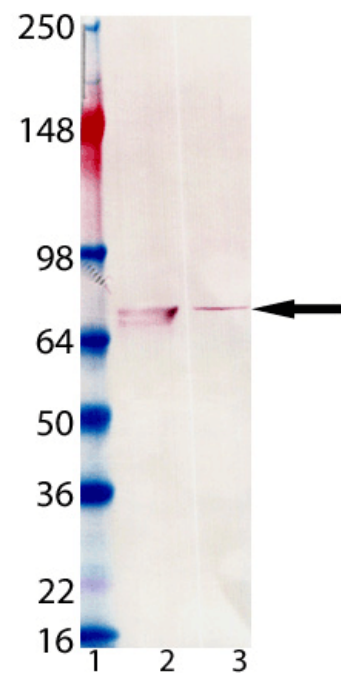
Immunohistochemistry analysis of human tonsil tissue using Zap-70 mAb (11H40), (Prod. No. ENZ-ABS709), at a dilution of 1:50. Panel A: Isotype Control, Panel B and C: ENZ-ABS709, diluted to 20µg/mL.

B

Immunohistochemistry analysis of human tonsil tissue using Zap-70 mAb (11H40), (Prod. No. ENZ-ABS709), at a dilution of 1:50. Panel A: Isotype Control, Panel B and C: ENZ-ABS709, diluted to 20µg/mL.

C

Immunohistochemistry analysis of human tonsil tissue using Zap-70 mAb (11H40), (Prod. No. ENZ-ABS709), at a dilution of 1:50. Panel A: Isotype Control, Panel B and C: ENZ-ABS709, diluted to 20µg/mL.



Western Blot analysis of Zap-70 mAb (11H40) (Prod. No. ENZ-ABS709): Lane 1: MW Marker, Lane 2: Jurkat lysate, Lane 3: Rat Thymus lysate.

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

ζ-chain associated protein kinase

IHC (PS), WB

Detects a band of ~70kDa by Western blot.

11H40

Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.

Mouse

Recombinant full length human Zap-70.

IgG1κ

Protein G affinity purified.

Immunohistochemistry (1:50)Western Blot (1:1,000,
ECL)Suggested dilutions/conditions may not be available
for all applications.Optimal conditions must be determined
individually for each application.

Purified from hybridoma tissue culture supernatant.

Human, Rat

P43403

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)