

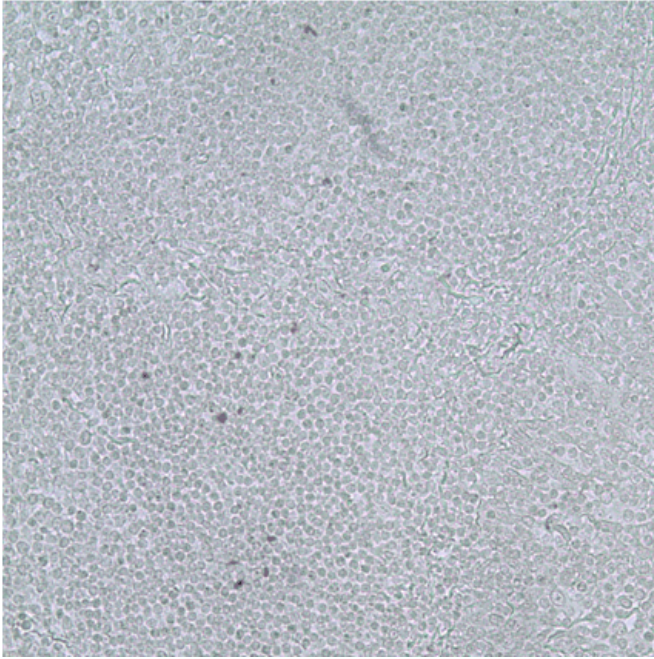
Zap-70 monoclonal antibody (11H40)

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.

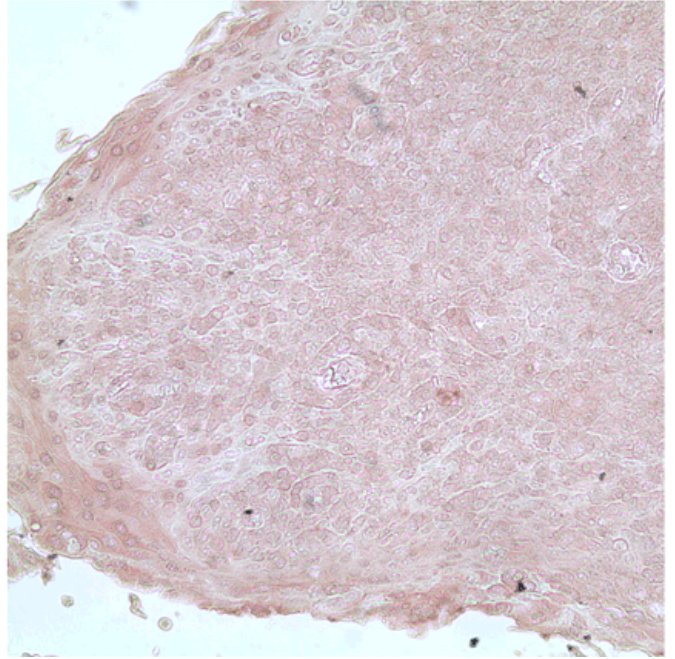
This antibody is covered by our [Worry-Free Guarantee](#).

Manuals, SDS & CofA

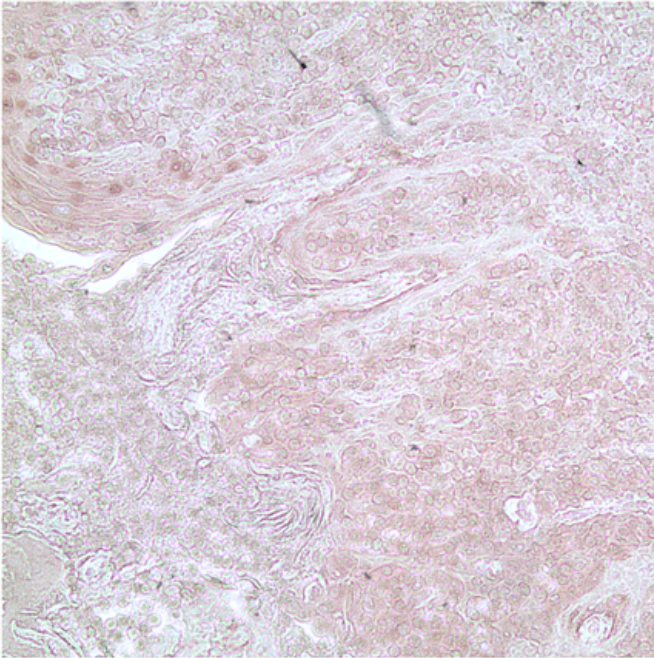
[View Online »](#)

A

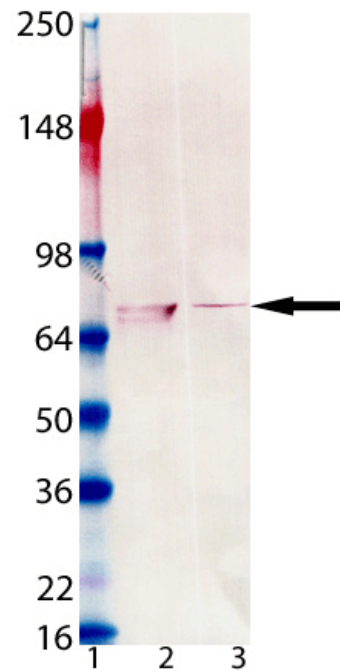
Immunohistochemistry analysis of human tonsil tissue using Zap-70 mAb (11H40), (Prod. No. ENZ-ABS709), at a dilution of 1:50. Panel A: Isotype Control, Panel B and C: ENZ-ABS709, diluted to 20µg/mL.

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 & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name ζ-chain associated protein kinase

Application IHC (PS), WB

Application Notes Detects a band of ~70kDa by Western blot.

Clone 11H40

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

Host Mouse

Immunogen Recombinant full length human Zap-70.

Isotype IgG1κ

Purity Detail Protein G affinity purified.

Recommendation Dilutions/Conditions 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.

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Human, Rat

UniProt ID P43403

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)