

c-Kit polyclonal antibody

c-Kit, also known as CD117 and Stem cell factor receptor, is a 145 kDa transmembrane tyrosine kinase encoded by the c-Kit protooncogene. c-Kit regulates a variety of biological responses including cell proliferation, apoptosis, chemotaxis, and adhesion. Ligand binding to the extracellular domain leads to autophosphorylation on several tyrosine residues within the cytoplasmic domain, and activation. c-Kit mutations correlate with tumor growth and progression in a variety of cancers including mast cell disease, gastrointestinal stromal tumor, acute myeloid leukemia, Ewing sarcoma, and lung cancer.

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

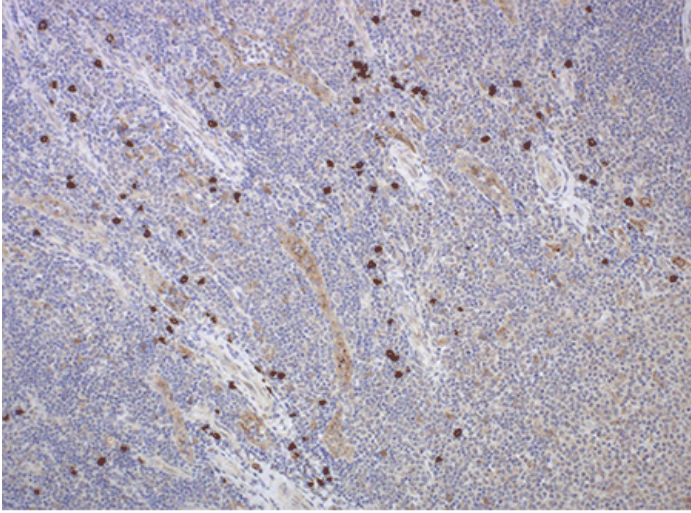
Ordering Information

[Order Online »](#)

ADI-KAP-TK005-E	100µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Formalin-fixed paraffin-embedded tonsil tissue stained using Enzo's C-KIT (CD117) antibody (Prod. No. ADI-KAP-TK005). IHC staining was performed on an automated Leica BOND III.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name CD117

Application IHC (PS), WB

Formulation Liquid. In PBS containing 50% glycerol.

Host Rabbit

Immunogen Synthetic peptide for C-terminal of human c-Kit.

Isotype IgG

Purity Detail Protein A-affinity purified.

Recommendation Dilutions/Conditions Immunohistochemistry (paraffin sections, 1:200)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.

Species Reactivity Human

UniProt ID P10721

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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)