

IL-17A polyclonal antibody

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

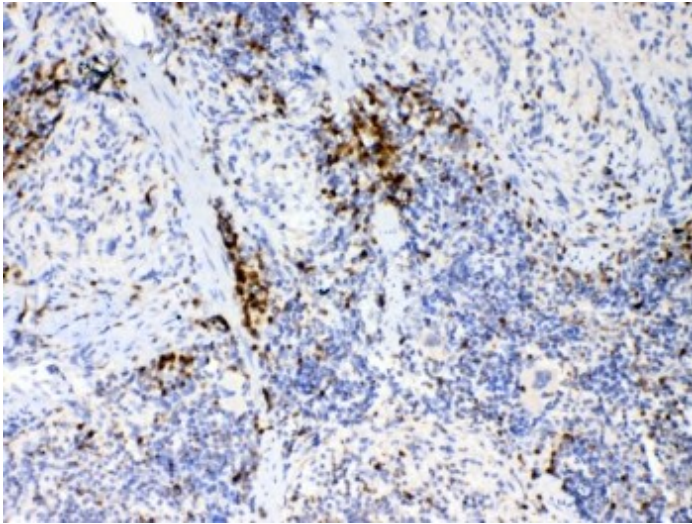
Ordering Information

[Order Online »](#)

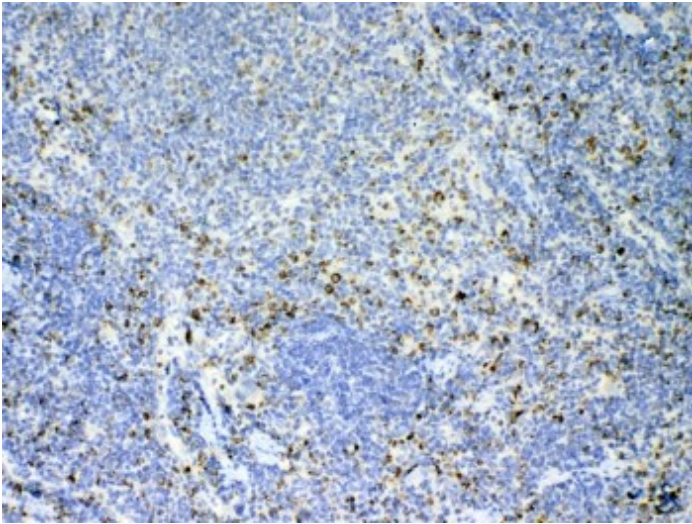
ENZ-ABS896-0100	100µg
-----------------	-------

Manuals, SDS & CofA

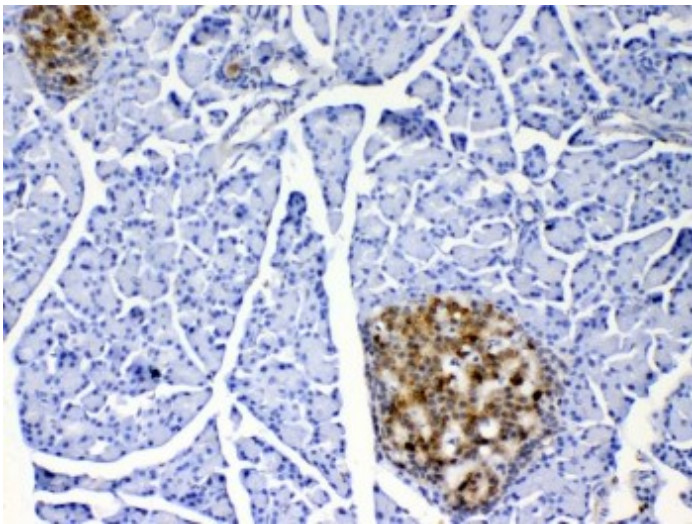
[View Online »](#)



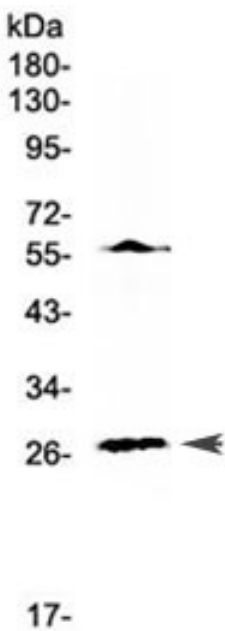
IHC testing of FFPE mouse spleen tissue with IL17A antibody at 1µg/ml. Required HIER: steam section in pH6 citrate buffer for 20 min and allow to cool prior to testing.



IHC testing of FFPE rat spleen tissue with IL17A antibody at 1µg/ml. Required HIER: steam section in pH6 citrate buffer for 20 min and allow to cool prior to testing.



IHC testing of FFPE rat pancreas tissue with IL17A antibody at 1µg/ml. Required HIER: steam section in pH6 citrate buffer for 20 min and allow to cool prior to testing.



Western blot testing of human Jurkat cell lysate with IL17A antibody at 0.5µg/ml. Expected molecular weight: 15-20 kDa (monomer), 30-40 kDa (homodimer), depending on glycosylation level.

Handling & Storage

Use/Stability	After reconstitution, the IL17A antibody can be stored for up to one month at +4°C. For long-term, aliquot and store at -20°C.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Interleukin-17A, Cytotoxic T-lymphocyte-associated antigen 8, CTLA-8
Application	ELISA, IHC (PS), WB
Formulation	Lyophilized from 1X PBS containing 2% Trehalose
Host	Rabbit
Immunogen	Recombinant mouse IL-17A (aa 26-158).
Isotype	IgG
Purity Detail	Antigen affinity purified.
Recommendation Dilutions/Conditions	ELISA (Capture) (0.1-0.5µg/ml)Immunohistochemistry (FFPE) (1-2µg/ml)Western blot (0.5-1µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse, Rat
UniProt ID	Q62386
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .

Last modified: July 12, 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)