

CD7 monoclonal antibody (CBC.37.8)

CD7 is a transmembrane glycoprotein expressed on T cells and NK cells. It plays a role in T-cell signaling and is used as a marker in leukemia and lymphoma diagnostics.

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

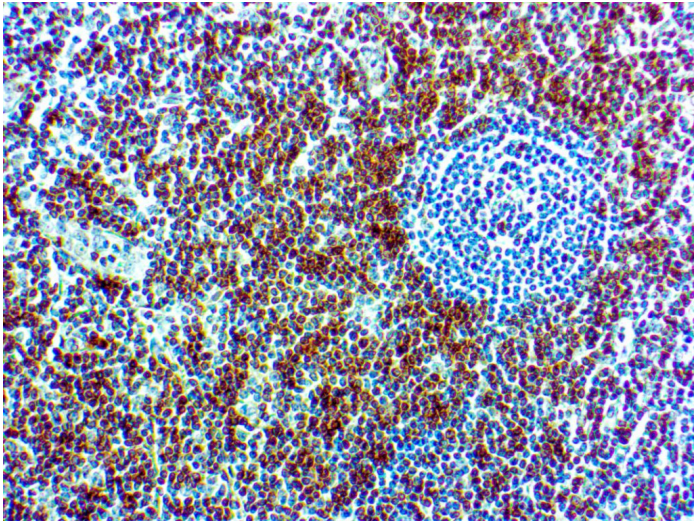
Ordering Information

[Order Online »](#)

ENZ-ABS1085-0500	500µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Formalin fixed paraffin embedded human tonsil stained with CD7 antibody

Handling	Avoid freeze/thaw cycles.
Long Term Storage	+4°C
Shipping	Blue Ice

Alternative Name	GP40
Application	IHC (PS)
Clone	CBC.37.8
Formulation	Contains sodium azide as a preservative.
Host	Mouse
Immunogen	A T lymphoblastoid cell line, established from a patient with acute lymphoblastic leukemia.
Isotype	IgG2bk
Positive Control	Tonsil
Purity Detail	Purified immunoglobulin.
Recommendation Dilutions/Conditions	Immunohistochemistry (paraffin sections) (1:50-1:200) Immunocytochemistry/Immunofluorescence (1:20-1:100) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.
Species Reactivity	Human
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .

Last modified: June 6, 2025



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)