

CD22 recombinant monoclonal antibody (340)

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

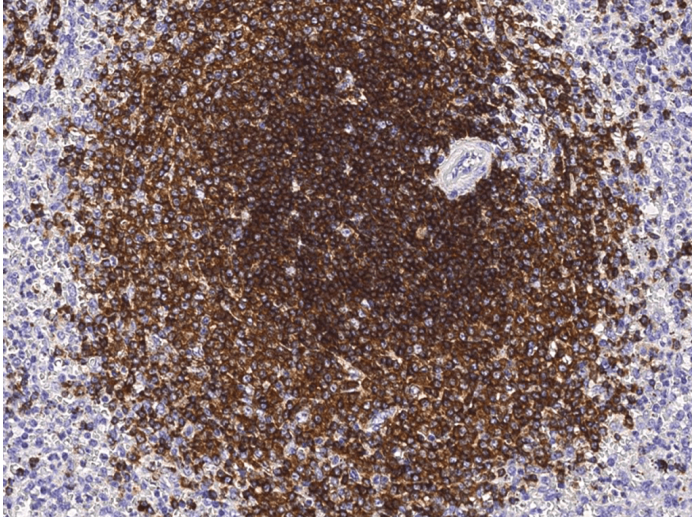
Ordering Information

[Order Online »](#)

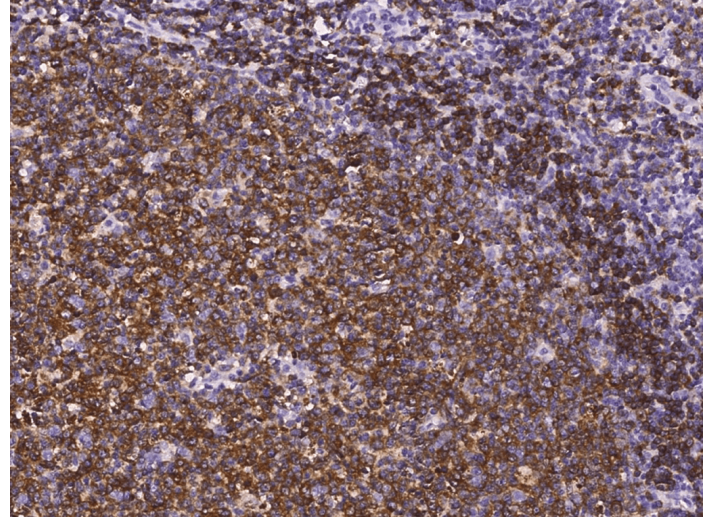
ENZ-ABS986-0100	100µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunochemical staining of human CD22 in human spleen with rabbit monoclonal antibody at 1:200 dilution, formalin-fixed paraffin embedded sections.



Immunochemical staining of human CD22 in human tonsil with rabbit monoclonal antibody at 1:200 dilution, formalin-fixed paraffin embedded sections.

Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	B-cell receptor CD22, B-lymphocyte cell adhesion molecule, BL-CAM, Sialic acid-binding Ig-like lectin 2, Siglec-2, T-cell surface antigen Leu-14
Application	ELISA, IHC (PS)
Clone	340
Formulation	Liquid. 0.2 µm filtered solution in PBS. Preservative free.
Host	Rabbit
Immunogen	Recombinant human CD22 protein.
Isotype	IgG
Purity Detail	Protein A purified.
Recommendation Dilutions/Conditions	ELISA (1:5'000-1:10'000)Immunohistochemistry (paraffin sections) (1:100-1:500)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



ENZO LIFE SCIENCES,
INC.
Phone: 800 942 0430
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(EUS) AG
Phone: +41 61 926 8989
info-
eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-
be@enzolifesciences.com

France
Phone: +33 472 440 655
info-
fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-
de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-
uk@enzolifesciences.com