

CD19 recombinant monoclonal antibody (109)

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

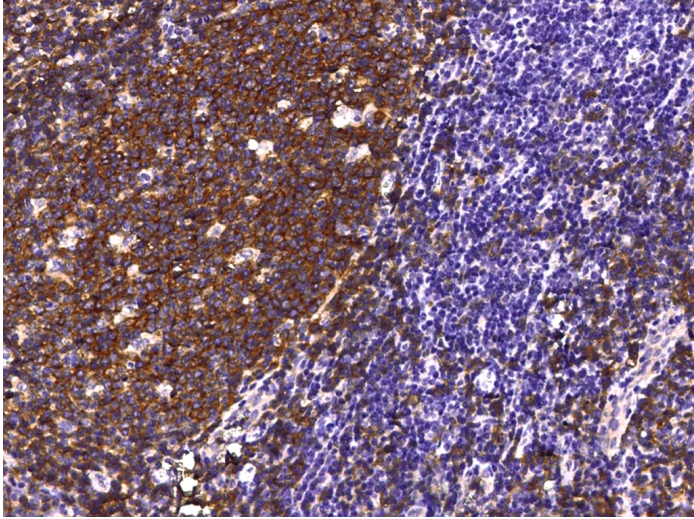
Ordering Information

[Order Online »](#)

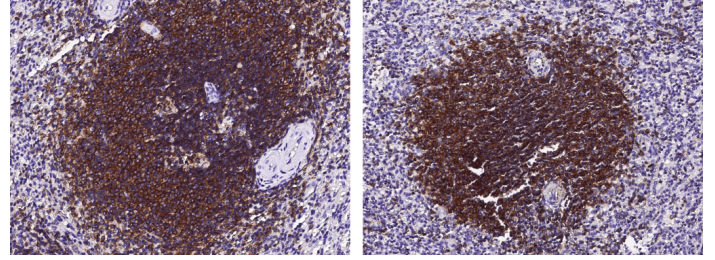
ENZ-ABS985-0100	100µl
-----------------	-------

Manuals, SDS & CofA

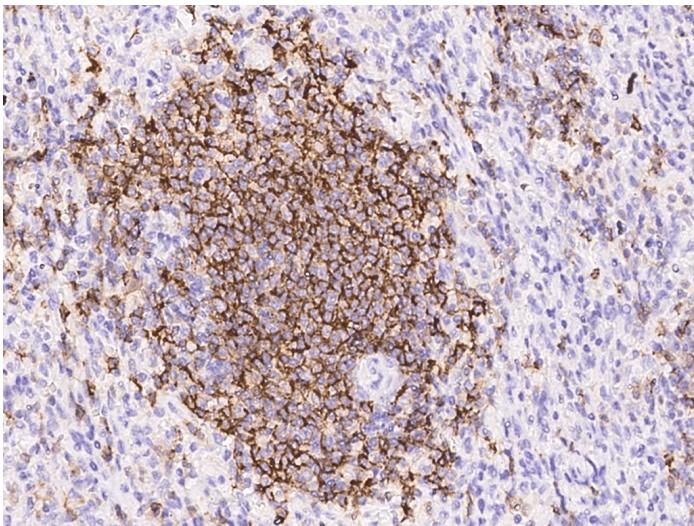
[View Online »](#)



Immunohistochemical staining CD19 in human tonsil with rabbit monoclonal antibody (1:300, formalin-fixed paraffin embedded sections). Positive staining was localized to membrane of B cells.



Immunohistochemical staining CD19 in human spleen with rabbit monoclonal antibody (1:300, formalin-fixed paraffin embedded sections). Positive staining was localized to membrane of B cells.



Immunohistochemical staining CD19 in human spleen with rabbit monoclonal antibody at 1:300 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-lymphocyte antigen CD19, B-lymphocyte surface antigen B4, Differentiation antigen CD19<T-cell surface antigen Leu-12

Application IHC (PS)

Clone 109

Formulation Liquid. 0.2 µm filtered solution in PBS. Preservative free.

Host Rabbit

Immunogen Recombinant human CD19 protein.

Isotype IgG

Purity Detail Protein A purified.

Recommendation Dilutions/Conditions 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](#)