CD22 recombinant monoclonal antibody (340)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

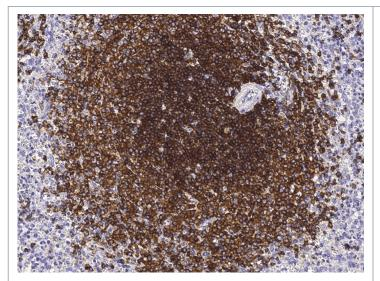
Order Online »

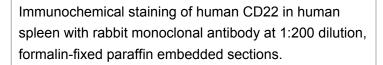
ENZ-ABS986-0100

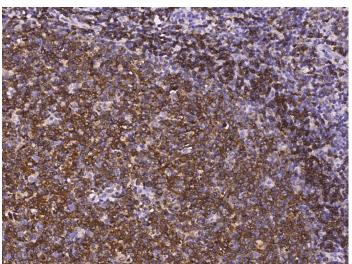
100µl

Manuals, SDS & CofA

View Online »







Immunochemical staining of human CD22 in human 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.

lgG lgG

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

.



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

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

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