CD32b recombinant monoclonal antibody (110)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

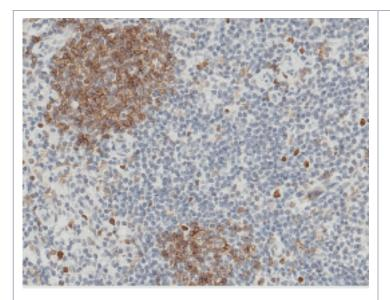
Order Online »

ENZ-ABS942-0100

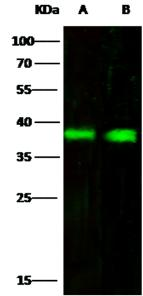
100µl

Manuals, SDS & CofA

View Online »



Immunohistochemical staining of formalin fixed, paraffinembedded human lymphonode showing membrane staining of B cells and some other mononuclear cell (1:10000).



Anti-CD32b rabbit monoclonal antibody at 1:500 dilution Lane A: Daudi Whole Cell Lysate Lane B: THP1 Whole Cell Lysate: Lysates/proteins at 30 µg per lane. Secondary: Goat Anti-Rabbit IgG H&L (Dylight800) at 1/10000 dilution. Developed using the Odyssey technique. Performed under reducing conditions. Observed band size:38 kDa

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 FCGR2B, Low affinity immunoglobulin gamma Fc region

receptor II-b

Application IHC (PS), WB

Clone 110

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

Host Rabbit

Immunogen Recombinant human CD32b protein.

lgG lgG

Purity Detail Protein A purified.

Recommendation Dilutions/ConditionsWestern Blot (1:500-1:1'000)Immunohistochemistry

(paraffin sections) (1:2500-1:15'000)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