CD79b recombinant monoclonal antibody (IGB/3170R)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

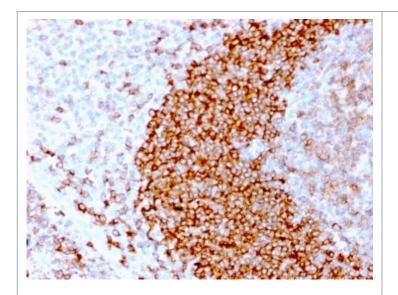
Order Online »

ENZ-ABS882-0100

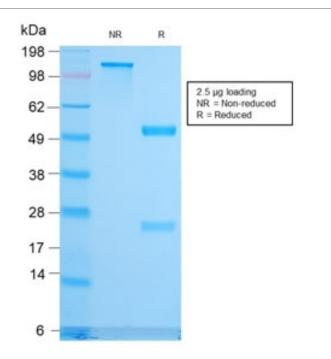
100µg

Manuals, SDS & CofA

View Online »



IHC staining of FFPE human tonsil tissue with recombinant CD79b antibody (clone IGB/3170R). HIER: boil tissue sections in pH 9 10mM Tris with 1mM EDTA for 10-20 min followed by cooling at RT for 20 min.



SDS-PAGE analysis of purified, BSA-free recombinant CD79b antibody (clone IGB/3170R) as confirmation of integrity and purity.

Handling & Storage

Handling Store the recombinant CD79b antibody at 2-8°C (with azide) or aliquot and store at -

20°C or colder (without azide).

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name B-cell antigen receptor complex-associated protein β

chain, Ig-β, B-cell-specific glycoprotein B29

Application ELISA, IHC (PS)

Clone IGB/3170R

Formulation Liquid. 0.2 mg/ml in 1X PBS containing 0.1 mg/ml BSA

(US sourced), 0.05% sodium azide.

Host Rabbit

Immunogen Recombinant human CD79b (aa 29-159).

Isotype IgGk

Purity Detail Protein A affinity chromatography purified.

Recommendation Dilutions/Conditions Immunohistochemistry (FFPE) (1-2µg/ml)Suggested

dilutions/conditions may not be available for all applications. Optimal conditions must be determined

individually for each application.

Species Reactivity Human

UniProt ID P40259

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

Last modified: July 12, 2024



info-

eu@enzolifesciences.com