

CD79b recombinant monoclonal antibody (IGB/3170R)

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

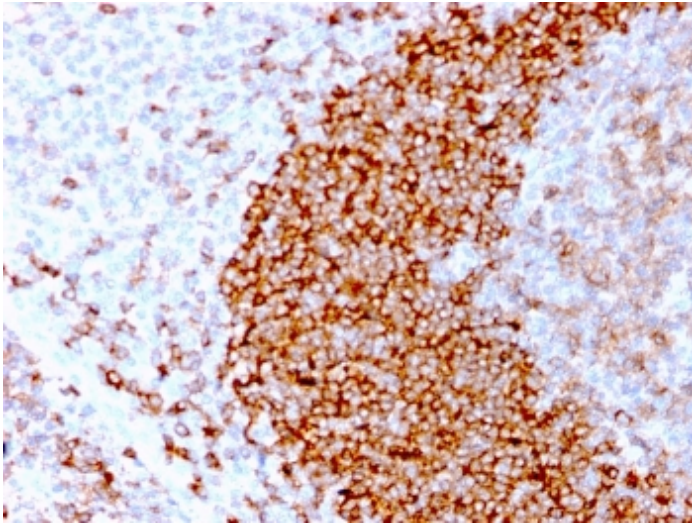
Ordering Information

[Order Online »](#)

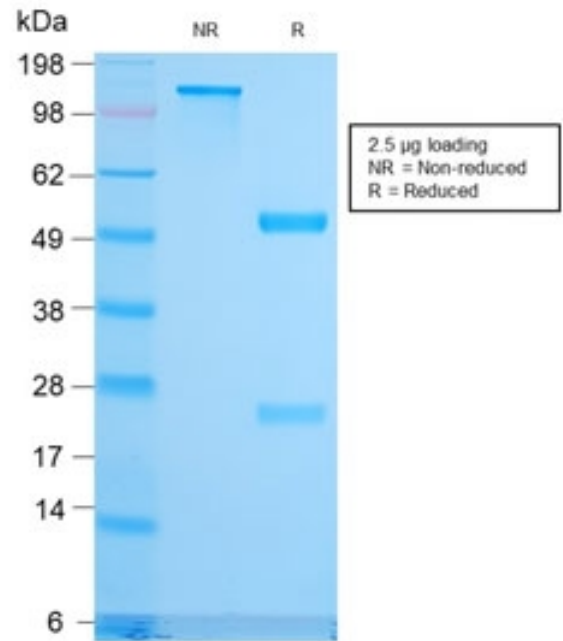
| | |
|-----------------|-------|
| ENZ-ABS882-0100 | 100µg |
|-----------------|-------|

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) (1-2 μ g/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](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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