

Bcl6 monoclonal antibody (SPM602)

Antibody to bcl-6 is helpful in a number of diagnostic settings: (1) In the differential diagnosis of small B-cell lymphoma. Follicular lymphoma will show bcl-6 (and CD10) positivity whereas other small B-cell lymphomas are usually negative. (2) Bcl-6 is an important prognostic marker in diffuse large B-cell lymphomas (DLBCL), where CD10, bcl-6 and MUM1/IRF4 are used to identify germinal center and activated B-cell phenotypes. (3) Bcl-6 can be valuable in distinguishing classical Hodgkin lymphoma from nodular lymphocyte predominant Hodgkin lymphoma (NLPHL). The Reed-Sternberg cells of classical Hodgkin lymphoma are bcl-6 negative whereas the large (L&H) cells of NLPHL are bcl-6 positive. In contrast, anti-Bcl-6 rarely stains mantle-cell lymphoma and MALT lymphoma.

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

Citations: 1

[View Online »](#)

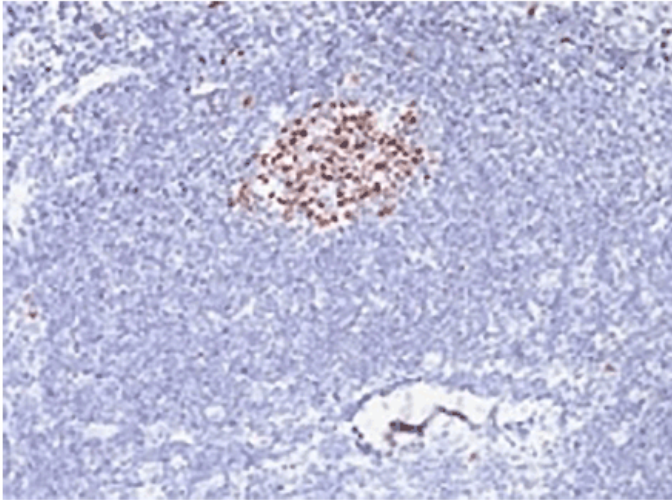
Ordering Information

[Order Online »](#)

ENZ-ABS771-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



IHC testing of formalin-fixed, paraffin-embedded human tonsil stained with anti-Bcl6 antibody

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name B-cell lymphoma 6 protein, B-cell lymphoma 5 protein, Bcl5

Application IHC

Clone SPM602

Formulation Liquid. In 1X PBS containing 0.1 mg/ml BSA and 0.05% sodium azide.

Host Mouse

Immunogen Recombinant human Bcl6.

Isotype IgG1κ

Purity Detail Protein G affinity purified.

Species Reactivity Human, Mouse

Specificity Recognizes a protein of 95kDa, which is identified as Bcl-6.

UniProt ID P41182

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

.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
Last modified: May 29, 2024
usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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