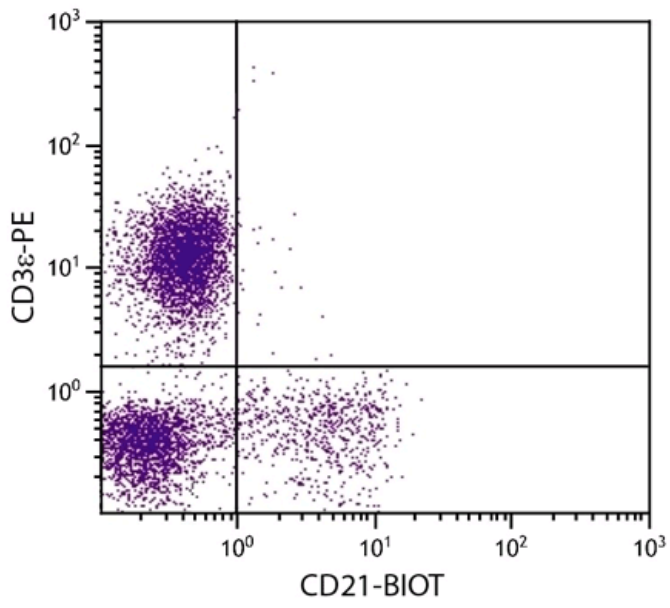


# CD21 monoclonal antibody (BB6-11C9.6) (Biotin)

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

Manuals, SDS & CofA

[View Online »](#)



Flow cytometry using CD21 Antibody [BB6-11C9.6]  
(Biotin)

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** +4°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Complement receptor 2, Cr2, Complement C3d receptor, Epstein-Barr virus receptor, EBV receptor

**Application** Flow Cytometry

**Clone** BB6-11C9.6

**Formulation** Liquid. In PBS containing sodium azide.

**Host** Mouse

**Isotype** IgG1

**Purity Detail** Purified.

**Species Reactivity** Porcine

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

**Last modified: May 29, 2024**



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-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

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