

# Complement component C4c monoclonal antibody (99-72-18)

C4c is a component in the Complement cascade and is a degradation product of C4b which is cleaved by C4b/C3b inactivator to yield C4d and C4c (soluble). C4c consists of a g-chain (30 kDa), a b-chain (75 kDa), and two a-domains denoted a3 (27 kDa) and a4 (14 kDa). C4c reacts with C1 and C2 to form C3 convertase of the classic activation pathway.

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

## Ordering Information

[Order Online »](#)

BPD-CAM-072-18-1	1mg
BPD-CAM-072-18-04	400µg

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Application ELISA, IF, IHC (PS), IP, WB

Clone 99-72-18

Formulation Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.

Host Mouse

Immunogen Synthetic peptide consisting of the last 15 C-terminal residues of the C4c a4-domain (Cys-EELVYELNPLDHRGR).

Isotype IgG2bk

Purity Detail Protein A/G-purified.

Source From culture supernatant.

Species Reactivity Human

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-  
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)