

# Complement component C5b-9 monoclonal antibody (aE11)

C5b-9, also known as the terminal complement complex (TCC), consists of C5b, C6, C7, C8 and C9, and it forms the membrane attack complex (MAC) as well as the non-lytic fluid-phase SC5b-9 complex (with protein S). The MAC forms channels in target cell membranes leading to cell lysis by osmotic leakage. The complexes contain neoantigens absent from the individual native components from which they are formed, and the antibody is directed against a neoepitope exposed on C9 when incorporated into the TCC.

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

Citations: 2

[View Online »](#)

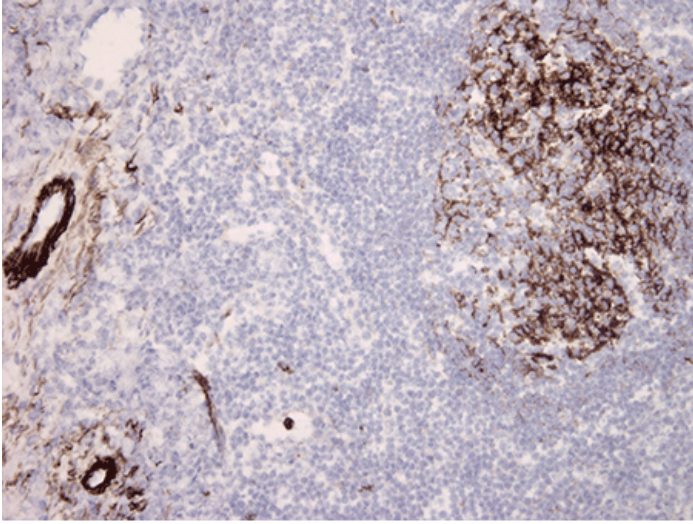
## Ordering Information

[Order Online »](#)

BPD-DIA-011-01-1	1mg
BPD-DIA-011-01-04	400µg

## Manuals, SDS & CofA

[View Online »](#)



IHC staining of frozen tonsil section. Vascular endothelia and dendritic cells of germinal center have strong and intensive label. Dilution: 1:10.

## Handling & Storage

Short Term Storage +4°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## Product Details

Alternative Name	Terminal complement complex, TCC
Application	ELISA, IF, IHC
Clone	aE11
Formulation	Liquid. In 10mM phosphate buffer, pH 7.4, containing 0.15M sodium chloride and 0.09% sodium azide.
Host	Mouse
Immunogen	Purified human complement component C5b-9.
Isotype	IgG2ak
Purity Detail	Protein A-affinity purified.
Species Reactivity	Baboon, Horse, Human, Porcine
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .

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-  
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)