

Complement component C5b-9 monoclonal antibody (aE11) (biotin conjugate)

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](#).

Ordering Information

[Order Online »](#)

BPD-DIA-011-01B-015

150µg

Manuals, SDS & CofA

[View Online »](#)

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

Clone aE11

Formulation Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.14M sodium chloride and 15mM sodium azide.

Host Mouse

Immunogen Purified human complement component C5b-9.

Isotype IgG2ak

Purity Detail Protein A/G-affinity purified.

Species Reactivity Baboon, Horse, Human, Porcine

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)