

β2-Microglobulin (caiman) monoclonal antibody (9-E10-C6)

Beta2-microglobulin is a 12-14kDa subunit of the major histocompatibility complex (MHC) class I. Beta2-microglobulin is a non-polymorphic molecule and is not involved in peptide binding, but is essential for MHC assembly. Beta2-microglobulin is also found in free form in the plasma. MHC class I is expressed on most nucleated cells in the caiman, including red blood cells.

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

Ordering Information

[Order Online »](#)

BPD-KSK-003-01-1	1mg
BPD-KSK-003-01-04	400µg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name β 2M

Application Affinity chromatography, Flow Cytometry, IP, WB

Clone 9-E10-C6

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

Host Mouse

Immunogen Recombinant caiman β 2-Microglobulin.

Isotype IgG1 κ

Purity Detail Protein A/G-purified.

Source From culture supernatant.

Species Reactivity Caiman

Specificity Reacts with free as well as MHC class I associated β 2M.

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-
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)