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 β2M

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 β2-Microglobulin.

lsotype lgG1κ

Purity Detail Protein A/G-purified.

Source From culture supernatant.

Species Reactivity Caiman

Specificity Reacts with free as well as MHC class I associated β2M.

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

Last modified: May 29, 2024

