
Product Information Sheet

Ordering Information

Product Number: V100010

Product Name: Anti-Beta-2-microglobulin (B2M) antibody *Mouse anti-human, monoclonal IgG1*

Unit Size: 50 ug

Storage Conditions: Freeze (<-15 °C)

Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Appearance: Solid

Molecular Weight: N/A

Soluble In: Water

Excitation Wavelength: N/A

Emission Wavelength: N/A

Application Notes

Human beta-2-microglobulin (B2M) is an 11.6 kDa non-glycosylated protein present in a variety of bodily fluids. It is a component of Class I MHC molecules and is necessary for the proper folding, binding of antigenic peptide, and surface expression of MHC-1 molecules. In long-term hemodialysis patients, B2M can accumulate to cause serious dialysis-related amyloidosis. Additionally, elevated levels of B2M are associated with rheumatoid arthritis, viral infections, cancer and aging. Anti-B2M monoclonal antibodies are highly specific for the detection of human B2M proteins. They can be used in a broad range of immunoassays such as latex enhanced turbidimetric immunoassays (LETIA), lateral-flow immunoassays (LFIA) and ELISA.