

Product Information Sheet

Ordering Information

Product Number: V100140

Product Name: Anti-Serum amyloid A protein antibody *Mouse anti-human, monoclonal IgG2b*

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

Serum amyloid A1 (SAA1) is a major acute-phase protein. It is produced primarily by hepatocytes in response to proinflammatory cytokines such as tumor necrosis factor α (TNF- α) or interleukin 6 (IL-6). Under inflammatory conditions, such as infection or malignancy, plasma concentrations of SAA1 can increase as high as 1000-fold. As such, it has become a promising reliable clinical indicator and biomarker for inflammatory diseases, chronic disorder and late-stage malignancy. Anti-SAA1 monoclonal antibodies are highly sensitive for the detection of human serum amyloid A1 immunogens. They can be used in the development of a broad range of immunoassays including gold immunochromatographic assays (GICA) and latex enhanced turbidimetric immunoassays (LETIA).