
Product Information Sheet

Ordering Information

Product Number: V100150

Product Name: Anti-C-reactive 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

C-reactive protein (CRP) is an annular, pentameric protein synthesized by the liver. It is present in blood plasma and is used primarily as a biomarker of inflammation. In response to inflammation, infection, or tissue damage, the level of CRP in human serum can increase 1000-fold. CRP also exhibits Ca²⁺ dependent binding to ligands, promote phagocytosis, and can induce production of hydrogen peroxide and inflammatory cytokines by monocytes. In clinical biology, CRP concentration can be used to evaluate the onset of cardiovascular disease. Anti-CRP monoclonal antibodies are highly specific for the detection of human C-reactive proteins. They can be used for a broad range of immunoassays such as lateral-flow immunoassay (LFIA), chemiluminescent immunoassay (CLIA), and latex enhanced turbidimetric immunoassays (LETIA).