

Product Information Sheet

Ordering Information

Product Number: V100135

Product Name: Anti-Interleukin-6 (IL6) 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

Interleukin 6 (IL-6) is a pleiotropic cytokine encoded by the IL6 gene in humans. It acts as both a pro-inflammatory cytokine and as an anti-inflammatory myokine. IL-6 is essential for stimulating acute phase protein synthesis, hybridoma growth, bone metabolism, inflammation and cancer progression. In almost all diseases involving chronic inflammation, such as diabetes rheumatoid arthritis and congestive heart failure, the concentration of IL-6 is elevated. Anti-IL-6 monoclonal antibodies are highly specific for the detection of human IL-6. They can be used for a broad range of immunoassays such as chemiluminescent immunoassays (CLIA) and rapid sandwich immunoassays.