
Product Information Sheet

Ordering Information

Product Number: V100030
Product Name: Anti-Cystatin-C protein 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

Cystatin C (CYSC) is a non-glycosylated, low molecular weight protein synthesized by most nucleated cells. It is present in all tissues and body fluids acting as a potent inhibitor of lysosomal proteinases and as an extracellular inhibitor of cysteine proteases. CYSC is primarily used as a biomarker of renal function, but also has associations with obesity, insulin resistance, hypertension, and cardiovascular mortality. Under normal conditions, CYSC in serum is 0.47 – 1.09 mg/L. When renal function is impaired, CYSC concentration in blood will change with glomerular filtration rate (GFR). During renal failure, GFR decreases while CYSC blood concentration increases more than 10-fold. Anti-CYSC monoclonal antibodies are highly sensitive for the detection of human CYSC proteins. They can be used for a broad range of CYSC immunoassays such as latex enhanced turbidimetric immunoassays (LETIA).