

Product Information Sheet

Ordering Information

Product Number:	V100170
Product Name:	Anti-Insulin-like growth factor-binding protein 1(IGFBP1) 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

Insulin-like growth factor-binding protein-1 (IGFBP-1), or placental protein 12 (PP12), belongs to a superfamily of IGFBPs that serve as carrier proteins for insulin-like growth factor (IGF-1). IGFBP-1 is synthesized and secreted by liver cells and by decidualized endometrial cells during pregnancy. It is expressed in liver, kidneys, and in amniotic fluid it is the most abundant IGFBP. IGFBP-1 binds equally to IGF-1 and IGF-II, and when phosphorylated IGFBP-1 exhibits higher binding affinities. In clinical biology, IGFBP-1 is a highly sensitive biomarker for the early diagnosis of premature rupture of fetal membranes (PROM) as the concentration of IGFBP-1 in amniotic fluid is 100 to 1000-fold higher than in human blood. Anti-IGFBP-1 monoclonal antibodies are highly sensitive for the detection of IGFBP-1 immunogens. They can be used for a broad range of immunoassays such as gold immunochromatography assays (GICA).