

Product Information Sheet

Ordering Information

Product Number: V100075

Product Name: Anti-Aspartate aminotransferase, cytoplasmic 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

Aspartate aminotransferase (AST) is a pyridoxal phosphate-dependent transaminase enzyme which exists in two isoenzymic forms mitochondrial (m-AST) and cytosolic (c-AST). It is essential for amino acid metabolism because it catalyzes the interconversion of aspartate and α -ketoglutarate to oxaloacetate and glutamate. AST is present in most organs with the highest concentrations found in the liver, heart, skeletal muscles, kidneys, brain and red blood cells. As a biomarker, AST has been used clinically to diagnose hepatitis, myocardial infarction, acute pancreatitis and musculoskeletal diseases. Anti-c-AST monoclonal antibodies are highly sensitive for the detection of human c-AST proteins. They can be used for a broad range of IHC applications and in immunoinhibition assay development.