

Product Information Sheet

Ordering Information

Product Number: V100110

Product Name: Anti-D-Dimer 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

D-dimer is a small protein fragment, or fibrin degradation product (FDP), present in blood after a blood clot has been degraded by fibrinolysis. It is composed of two D fragments of the fibrin protein joined together by a cross-link. Elevated d-dimer levels are a key indicator of thrombotic events, indicating excess fibrinolysis, following activation of coagulation. In clinical biology, it is used in the diagnosis of the blood disorder disseminated intravascular coagulation. Anti-D-Dimer monoclonal antibodies are highly sensitive for the detection of human D-dimer immunogens. They can be used in a broad range of immunoassays such as lateral-flow immunoassays (LFIA) and in IVD assay development.