
Product Information Sheet

Ordering Information

Product Number: V100130
Product Name: Anti-Trypsinogen-2 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

Trypsinogen-2 (PRSS2) is a zymogen of the pancreatic proteinase trypsin. It is synthesized by the pancreas and is present at high concentrations in the pancreatic juice. Trypsinogen-2 is secreted into the duodenal lumen of the small intestine, where it is activated by enterokinases to produce the protein digestive enzyme trypsin. In clinical biology, measurement of serum trypsinogen-2 is a useful tool in diagnosing acute pancreatitis and cystic fibrosis. Patients with such disorders express high levels of trypsinogen-2. Anti-PRSS2 monoclonal antibodies are highly-specific for the detection of human trypsinogen-2 immunogens. They can be used for a broad range of immunoassays such as ELISA, as well as in the development of IVD immunoassays.