

Anti-Procalcitonin (PCT) antibody *Mouse anti-human, monoclonal IgG2b*

Catalog Number: V100115, V100116

Unit Size: 50 ug, 1 mg

Product Details

Storage Conditions	Freeze (< -15 °C)
Expiration Date	12 months upon receiving

Biological Properties

Appearance	Liquid
Soluble In	Water

Applications

Procalcitonin (PCT) is a peptide precursor of calcitonin and is composed of 116 amino acids. It is produced by the C cells of the thyroid and by the neuroendocrine cells of the lung and the intestine. Under normal conditions, the level of PCT in the blood of healthy individuals is very low (< .05 ng/mL) with a half-life of 25 to 30 hours in serum. In response to a pro-inflammatory stimulus, such as a pathogen, the level of PCT may increase as high as 100 µg/mL making it a suitable biomarker of bacterial infection. In clinical biology, PCT has shown to be a useful tool in the diagnosis of sepsis. Anti-PCT monoclonal antibodies are highly specific for the detection of human procalcitonin. They can be used for a broad range of immunoassays such as chemiluminescent immunoassays (CLIA).