

Product Information Sheet

Ordering Information

Product Number: V100115

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

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

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 ?g/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).