

Product Information Sheet

Ordering Information

Product Number: V100000

Product Name: Anti-Neutrophil gelatinase-associated lipocalin (NGAL) 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

Neutrophil gelatinase-associated lipocalin (NGAL) is a protein expressed mostly in neutrophils. It is also present in lower levels in the kidney, prostate and epithelia of the respiratory and alimentary tracts. NGAL is a vital component in the innate immune response to bacterial infection. It limits bacterial growth by sequestering iron-containing microorganisms. NGAL's small size and protease resistant properties makes it easily excreted and detected in the urine. Under normal conditions, NGAL levels in the urine and plasma are low. However, in response to kidney injury NGAL levels increase sharply reaching diagnostic levels quickly. In clinical biology, NGAL is a novel biomarker of acute kidney injury (AKI) and is used in the early diagnosis of chronic kidney disease. Anti-NGAL monoclonal antibodies are highly sensitive for the detection of human NGAL. They can be used for a broad range of immunoassays such as chemiluminescent immunoassays (CLIA), latex enhanced turbidimetric immunoassays (LETIA), lateral-flow immunoassays (LFIA) and ELISA.