

Product Information Sheet

Ordering Information

Product Number: V100165

Product Name: Anti-Prolactin (PRL) 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

Prolactin (PRL), also known as luteotropic hormone, is a protein secreted from the pituitary gland. It plays a vital role in metabolism, immune system regulation and pancreatic development, but is best known for its role in producing milk in mammals. In normal individuals, PRL concentrations increase in response to physiologic stimuli such as eating, mating, exercise, stress and hypoglycemia. PRL levels are also elevated during pregnancy, lactation, postpartum and in newborn infants. In clinical biology, prolactin tests are used in the diagnosis of prolactinoma, a prolactin-producing tumor of the pituitary gland. Anti-PRL monoclonal antibodies are highly sensitive for the detection of human purified prolactin. They can be used for a broad range of immunoassays such as chemiluminescent immunoassays (CLIA) and in the development of highly sensitive sandwich immunoassays.