

Product Information Sheet

Ordering Information

Product Number: V100180

Product Name: Anti-Follicle-stimulating hormone 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

Follicle-stimulating hormone (FSH) is a 35.5 kDa glycoprotein polypeptide hormone synthesized and secreted by the gonadotropic cells of the anterior pituitary gland. It is essential for regulating the development, growth, pubertal maturation and reproductive processes of the body. In clinical biology, FSH levels can be used to evaluate physiological disorders. For instance, FSH and luteinizing hormone levels are elevated in primary gonadal failure, primary hypogonadism, complete testicular feminization syndromes and menopause. In pituitary or hypothalamus FSH and LH levels are both decreased. Anti-FSH monoclonal antibodies are highly sensitive for the detection of human follicle-stimulating hormones. They can be used for a broad range of immunoassays such as lateral-flow immunoassays (LFIA), chemiluminescence immunoassays (CLIA) and rapid sandwich immunoassays.