
Product Information Sheet

Ordering Information

Product Number: V100060

Product Name: Anti-Epididymal secretory protein 4 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

Human epididymis secretory protein E4 (HE4) is a protein encoded by the WFDC2 gene in humans. This gene is expressed primarily in pulmonary epithelial cells, but has found to be expressed in some ovarian cancers. As such, HE4 serves as a serum biomarker in women with a pelvic mass to distinguish benign from malignant disease and as a tumor marker for ovarian cancer. Anti-HE4 monoclonal antibodies are highly sensitive for the detection of human HE4. They can be used for a broad range of immunoassays such as chemiluminescence immunoassay (CLIA) and rapid sandwich immunoassays.