

## Product Information Sheet

### **Ordering Information**

Product Number: V100220

Product Name: Anti-Cortisol 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**

Cortisol is a steroid hormone of the glucocorticoid class that is synthesized from cholesterol by the adrenal gland. While it has a multiple of roles, Cortisol is best known for its effects on metabolism, its immune system function, its ability to increase blood sugar via gluconeogenesis and as a precursor for other steroid hormones. Cortisol levels are regulated by the hypothalamic-pituitary-adrenal complex. In response to the release of corticotropin-releasing hormone (CRH) by the hypothalamus, adjacent cells of the anterior pituitary secrete an adrenocorticotrophic hormone (ACTH) into the vascular system. Blood transports ACTH to the adrenal cortex where it stimulates the synthesis of cortisol. In clinical biology, measuring cortisol levels is a useful biomarker for Cushing's disease, which is characterized by the overproduction of cortisol. Additionally, abnormal cortisol levels have also been implicated in prostate cancer, depression and schizophrenia. Anti-cortisol monoclonal antibodies are highly sensitive for the detection of cortisol immunogens. They can be used for a broad range of immunoassays such as ELISA