

**Anti-Estradiol antibody \*Mouse anti-human, monoclonal IgG1\***Catalog number: V100185, V100186  
Unit size: 50 ug, 1 mg**Product Details**

---

Storage Conditions	Freeze (<-15 °C)
Expiration Date	12 months upon receiving

**Chemical Properties**

---

Appearance	Solid
Soluble In	Water

**Applications**

---

Estradiol (E2) is an estrogen steroid hormone produced primarily within the follicles of ovaries, but is also produced in the testicles, adrenal glands, breasts and brain. It is the major female sex hormone responsible for the regulation of estrous and menstrual reproductive cycles, and is necessary for female development. It is also produced in males, at considerably lower levels compared to females, as an active metabolic product of testosterone. In clinical biology, measuring serum estradiol levels is used to assess the activity of female ovaries, to detect baseline estrogen in females with menstrual dysfunction, and to detect menopause. Anti-estradiol monoclonal antibodies are highly sensitive for the detection of estradiol immunogens. They can be used for a broad range of immunoassays such as ELISA.