

**Anti-17-alpha-hydroxyprogesterone
antibody *Mouse anti-human, monoclonal
IgG1***Catalog number: V100195, V100196
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

17 alpha-hydroxyprogesterone (17 α -OHP) is a steroid hormone synthesized from progesterone. It acts as an agonist of the progesterone receptor and is a precursor in the biosynthesis of many other steroids such as glucocorticoids, androgens and estrogens. In clinical biology, measuring 17 α -OHP levels in the blood are useful in the detection and evaluation of congenital adrenal hyperplasia (CAH). Patients with CAH express insufficient amounts of the enzyme 21-hydroxylase, which is necessary for the conversion of 17 α -OHP into cortisol. As a result, 17 α -OHP levels in the blood build up. Anti-17 α -OHP monoclonal antibodies are highly sensitive for the detection of 17 α -OHP. They can be used for a broad range of immunoassays such as competitive chemiluminescence immunoassays (CLIA) and ELISA.