

**Purified Mouse Anti-human  $\beta$ -hCG Antibody**  
**\*HCG-61, monoclonal\***Catalog number: V103225  
Unit size: 0.1 mg**Product Details**

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	Lot specific (please consult certificate of analysis for given lot)
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 15 mM sodium azide, 0.2% (w/v) BSA

**Antibody Properties**

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	$\beta$ -hCG
Clone	HCG-61

**Biological Properties**

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	ICC, ELISA

**Applications**

Choriogonadotropin subunit  $\beta$  3 (also known as CG- $\beta$ , Choriogonadotropin subunit  $\beta$  or Chorionic gonadotropin chain beta) is a protein with a molecular weight of 21 kDa, expressed in the cytoplasm, extracellular region and extracellular space of cells. In humans, choriogonadotropin subunit  $\beta$  3 has been closely linked to vital functions such as hormone activity. It is the subject of comprehensive application in part because of the fact that it acts in the G protein-coupled receptor signaling pathway and hormone-mediated signaling pathway. It is thought to be essential to transcription by RNA polymerase II. Choriogonadotropin subunit  $\beta$  3 has been found to be involved in organismal processes, in particular, cell-cell signaling, apoptotic process and peptide hormone processing.