

**Biotin Mouse Anti-human C-peptide
Antibody *C-PEP-01, monoclonal***Catalog number: V103370
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	C-peptide
Clone	C-PEP-01
Conjugate	Biotin

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with Biotin under optimal conditions
Application	ELISA, RIA, IHC(P), ICC

Applications

Proinsulin C-peptide (C-peptide) has shown evidence of use in a multitude of applications, particularly in the field of pharmacological research. It is vital to the examination of key biological diseases like autoimmune disease and dual-hormone-deficiency disease, specifically in cell-lines like vascular smooth muscle cells, assess pancreatic beta cells and L6 myoblasts. Additionally, it has been widely used in the study of vital biological activities like ATPase activity, HGF activity and Metal-Modulated Activity. It has a variety of applications in the study of important cell-based enzymes like Akt kinase, normal co-release and serine protease.