

**Anti-Fatty acid-binding protein antibody**  
**\*Mouse anti-human, monoclonal IgG1\***

Catalog number: V100105, V100106

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**

---

Heart-type fatty acid-binding protein (H-FABP) is a 15 kDa cytoplasmic protein encoded by the FABP3 gene in humans. It is involved in fatty acid metabolism transporting fatty acids from the cell membrane to the mitochondria for oxidation. Following ischemic episodes, H-FABP is released from cardiac myocytes making it a sensitive biomarker for myocardial infarction. H-FABP is 20 times more specific to cardiac muscle than myoglobin and is found 10-fold lower levels in skeletal muscle than heart muscle. Anti-H-FABP monoclonal antibodies are highly sensitive for the detection of human H-FABP. They can be used for a broad range of immunoassays such as lateral-flow immunoassays (LFIA), chemiluminescent assays (CLIA) and sandwich immunoassays.