

Purified Mouse Anti-human MFG Antibody
MFG-06, monoclonalCatalog number: V1031750
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	MFG
Clone	MFG-06

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	WB, IHC(P)

Applications

Lactadherin is a 50 kDa protein that can be found in the extracellular space, endoplasmic reticulum lumen and membrane of cells. It is also known as HMFG, breast epithelial antigen BA46 and MFGM. In humans, MFGM takes part in processes such as organismal processes, for example, phagocytosis, recognition, post-translational protein modification and phagocytosis, engulfment, and moreover, binds to phosphatidylethanolamine, integrin and phosphatidylserine. Sequencing of MFGM has supported it contains 3 conserved structural units: F5/8 type C 2, EGF-like and F5/8 type C 1 domain. It is an enhancer of phagocytosis.