

**Purified Mouse Anti-human SUSD2
Antibody *W5C5, monoclonal***

Catalog number: V1032180

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	SUSD2
Clone	W5C5

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	ICC, FC (QC TESTED)

Applications

Sushi domain-containing protein 2 is a 90 kDa transmembrane protein that can be located in the integral component of membrane, plasma membrane and extracellular exosome of cells. Sequencing of Sushi domain-containing protein 2 has shown it contains 6 types of conserved structural units: VWFD, SMB, cytoplasmic, extracellular, AMOP and Sushi domain. Sushi domain-containing protein 2 inhibits cell cycle G1/S phase transition and cell division.