

**APC Mouse Anti-human SUSD2 Antibody**  
**\*W5C5, monoclonal\***

Catalog number: V1032185

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
Conjugate	APC

**Biological Properties**

---

Preparation	Antibody purified by affinity chromatography and then conjugated with APC under optimal conditions
Application	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.