

**Purified Mouse Anti-human AGR3 Antibody**  
**\*AGR3.2, monoclonal\***

Catalog number: V103055

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	AGR3
Clone	AGR3.2

**Biological Properties**

---

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

**Applications**

---

Anterior gradient protein 3 (also known as Protein disulfide isomerase family A, member 18, Breast cancer membrane protein 11 or Anterior gradient 3 homolog) is a protein with a molecular weight of 22 kDa, found in the intracellular membrane-bounded organelle and endoplasmic reticulum of cells. In Homo sapiens, anterior gradient protein 3 reacts with dystroglycan. It inhibits cell death.