

Purified Mouse Anti-human AGR3 Antibody
AGR3.1, monoclonalCatalog number: V103050
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	AGR3.1

Biological Properties

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