

**Purified Mouse Anti-human AGR2/AGR3
Antibody *AGR3.4, monoclonal***Catalog number: V103045
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	AGR2/AGR3
Clone	AGR3.4

Biological Properties

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

Applications

Anterior gradient protein 2 homolog (sometimes referred to as Secreted cement gland protein XAG-2 homolog or HPC8) is a protein with a molecular weight of 23 kDa, located in the extracellular space and endoplasmic reticulum of cells. In humans, AG-2 is thought to be essential to organismal processes, for example, lung goblet cell differentiation, mucus secretion and digestive tract morphogenesis. It binds with dystroglycan, identical protein and epidermal growth factor receptor. AG-2 is the subject of intensive examination because of the fact that it is a component of the positive regulation of epidermal growth factor receptor signaling pathway. Additionally, it acts to positively regulate protein localization to plasma membrane, gene expression and cell-substrate adhesion while also inhibits cell death.