

**Purified Mouse Anti-human/pig GFAP  
Antibody \*GF-02, monoclonal\***Catalog number: V103725  
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	Tris buffered saline (TBS) solution, 15 mM sodium azide, 0.2% (w/v) BSA

**Antibody Properties**

Species Reactivity	Human, pig
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	GFAP
Clone	GF-02

**Biological Properties**

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

**Applications**

Glial fibrillary acidic protein is a protein with a molecular weight of 50 kDa, found in the cytoplasm, cytosol and intermediate filament of cells. Sequencing of glial fibrillary acidic protein has exemplified it contains a primary structural unit, the IF rod domain. Glial fibrillary acidic protein aids in protein-containing complex assembly and chaperone-mediated autophagy. It reacts with identical protein. It plays an important role in organismal processes, for example, intermediate filament organization. Mutations and abnormalities in glial fibrillary acidic protein have been associated with a number of diseases, namely, Alexander disease (ALXDRD). Alexander disease, an autosomal dominant inheritance disorder characterized by hydrocephalus, progressive macrocephaly and increased CSF protein, has in particular been of interest to scientists.