

**FITC Mouse Anti-human/pig GAPDHS  
Antibody \*Hs-8, monoclonal\***Catalog number: V103715  
Unit size: 50 tests**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	Stabilizing 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	GAPDHS
Clone	Hs-8
Conjugate	FITC

**Biological Properties**

Preparation	Antibody purified by affinity chromatography and then conjugated with FITC under optimal conditions
Application	FC (QC TESTED)

**Applications**

Glyceraldehyde-3-phosphate dehydrogenase, testis-specific is a 45 kDa protein that can be found in the nucleus and cytosol of cells. It is alternatively called Spermatogenic glyceraldehyde-3-phosphate dehydrogenase and Spermatogenic cell-specific glyceraldehyde 3-phosphate dehydrogenase 2. In Homo sapiens, spermatogenic cell-specific glyceraldehyde 3-phosphate dehydrogenase 2 has been associated with critical functions such as glyceraldehyde-3-phosphate dehydrogenase (NAD<sup>+</sup>) (phosphorylating) activity, and also, plays a role in the upregulation of glycolytic process. Spermatogenic cell-specific glyceraldehyde 3-phosphate dehydrogenase 2 recognizes NADP and NAD, and also, takes part in organismal processes, in particular, glycolytic process and gluconeogenesis.