

**Purified Mouse Anti-human/pig GAPDHS
Antibody *Hs-8, monoclonal***Catalog number: V103710
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 | GAPDHS |
| Clone | Hs-8 |

Biological Properties

| | |
|-------------|--|
| Preparation | Antibody purified by affinity chromatography and then conjugated with under optimal conditions |
| Application | WB, ICC, 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⁺) (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.