

**Purified Mouse Anti-FITC Antibody**  
**\*Monoclonal\***Catalog number: V102105  
Unit size: 1.0 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	, 0.2% (w/v) BSA

**Antibody Properties**

Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	FITC

**Biological Properties**

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	Flow Cytometry (FACS)

**Applications**

Fluorescein is a fluorescent indicator. It has been widely used in the research of important cellular enzymes like protein kinase, dilute aqueous bases and DNA apurinic or apyrimidinic site lyase. Fluorescein has a variety of applications in the study of vital biological receptors like peptide ligand-binding receptors and thyrotropin releasing hormone receptors. Direct photolysis, continuous electrophoresis and harmonised basis are, for instance, just a few of the processes that it has helped researchers to better understand. It is essential for the research of critical biological activities like administrative activities. Moreover, it has a variety of applications in the examination of key biological proteins such as membrane bound protein and bloom syndrome protein.