

Purified Rabbit Anti-Calmodulin Binding Peptide Antibody *Polyclonal*

Catalog number: V102035

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	, 0.2% (w/v) BSA

Antibody Properties

Class	Primary
Clonality	Polyclonal
Host	Rabbit
Immunogen	Calmodulin Binding Peptide

Biological Properties

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

Applications

A 5.3 kDa peptide that binds to calmodulin (CaM) with high affinity, derived from smooth muscle myosin light-chain kinase (MLCK).