

HRP Rabbit Anti-Calmodulin Binding Peptide Antibody

Polyclonal

Catalog Number: V102050

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
Formulation	HEPES – saline with 0.002% Thimerosal and Stabilizer

Antibody Properties

Class	Primary
Clonality	Polyclonal
Host	Rabbit
Immunogen	Calmodulin Binding Peptide

Biological Properties

Appearance	Liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with HRP under optimal conditions
Application	FACS

Spectral Properties

Conjugate	HRP
-----------	-----

Applications

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