

**Purified Mouse Anti-human/mouse/rat
STIM1 Antibody *CDN3H4, monoclonal***Catalog number: V1032160
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	Phosphate-buffered saline (PBS, pH 7.2), 15 mM sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human, mouse, rat
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	STIM1
Clone	CDN3H4

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	IP, WB, IHC(P), ICC

Applications

This is an antibody that targets STIM1.