

**Purified Rabbit Anti-mouse/rat GABA(B)
R1a Antibody *PAb (549), monoclonal***Catalog number: V103685
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	Mouse, rat
Class	Primary
Clonality	Monoclonal
Host	Rabbit
Immunogen	GABA(B) R1a
Clone	PAb (549)

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	WB, ICC

Applications

This is an antibody that targets GABA(B) R1a.