Glycine receptor polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

BML-SA510-0200

200µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application IHC, WB

Formulation Lyophilized from 5mM ammonium bicarbonate.

Host Rabbit

Immunogen Synthetic peptide corresponding to a portion of N-terminal

rat glycine receptor subunit α1.

Species Reactivity Human, Mouse, Rat

UniProt ID P07727

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024

