P2Y14 receptor (extracellular) polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

BML-SA590-0050

50µl

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name P2RY14

Application Flow Cytometry, WB

Formulation Lyophilized.

Host Rabbit

Immunogen Synthetic peptide corresponding to aa 172-188 of human

P2RY14.

Species Reactivity Human, Mouse

UniProt ID Q15391

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

•

Last modified: May 29, 2024

