VE-cadherin polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Citations: 2

View Online »

Ordering Information

Order Online »

ENZ-ABS661-0100

100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Vascular endothelial cadherin, Cadherin-5, CD144

Application ICC, IF, IHC, IP, WB

Formulation Liquid. In 0.01M PBS, pH 7.4, containing 0.05% Proclin-

300 and 50% glycerol.

Host Rabbit

Immunogen Recombinant human VE-cadherin (aa 68-307).

Purity Detail Purified by antigen-specific affinity chromatography

followed by protein A affinity chromatography.

Species Reactivity Human

UniProt ID P33151

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

•

Last modified: May 15, 2025

