Filaggrin polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Citations: 3

View Online »

Ordering Information

Order Online »

ENZ-ABS181-0100

100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application IF, IHC, WB

Formulation Phosphate-buffered solution, pH 7.2, containing 0.09%

sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to aa 24-39 in the partial

sequence of filaggrin.

Purity Detail Affinity purified.

Species Reactivity Human, Mouse, Primate

Technical Info / Product Notes Cited samples:

For an overview on cited samples please click here.

UniProt ID P20930

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

.

Last modified: September 24, 2025

