

# NEDD8 polyclonal antibody

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

ENZ-ABS376-0025	25µl
-----------------	------

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** Neddylin, Neural precursor cell expressed developmentally down-regulated protein 8, Ubiquitin-like protein Nedd8

**Application** IHC, WB

**Formulation** Liquid. In PBS, pH 7.3, containing 0.05% sodium azide and 50% glycerol.

**Host** Rabbit

**Immunogen** Recombinant human NEDD8.

**Isotype** IgG

**Purity Detail** Affinity-purified.

**Recommendation Dilutions/Conditions** Immunohistochemistry (1:50-1:100)Western Blot (1:500-1:2,000)Optimal conditions must be determined individually for each application.

**Species Reactivity** Human, Mouse, Rat

**UniProt ID** Q15843

**Worry-free Guarantee** This antibody is covered by our [Worry-Free Guarantee](#)

.