

RAB5 polyclonal antibody

The Rab family of GTP-binding proteins are members of the Ras superfamily of small G proteins, and participate in multiple steps of the membrane trafficking. Rab3A participates in the fusion of synaptic vesicles during neurotransmitter release, a process regulated by GDP/GDP exchange cofactors such as GAP, GDI, GRF, and rabphilin. Rab4 regulates recycling of proteins from the early endosome to the surface of many cell types, regulating glucose transporter and transferrin receptor cell surface localization. Rab5 regulates the fusion of plasma membrane-derived clathrin-coated vesicles with early endosomes and homotypic fusion among early endosomes.

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

Citations: 4

[View Online »](#)

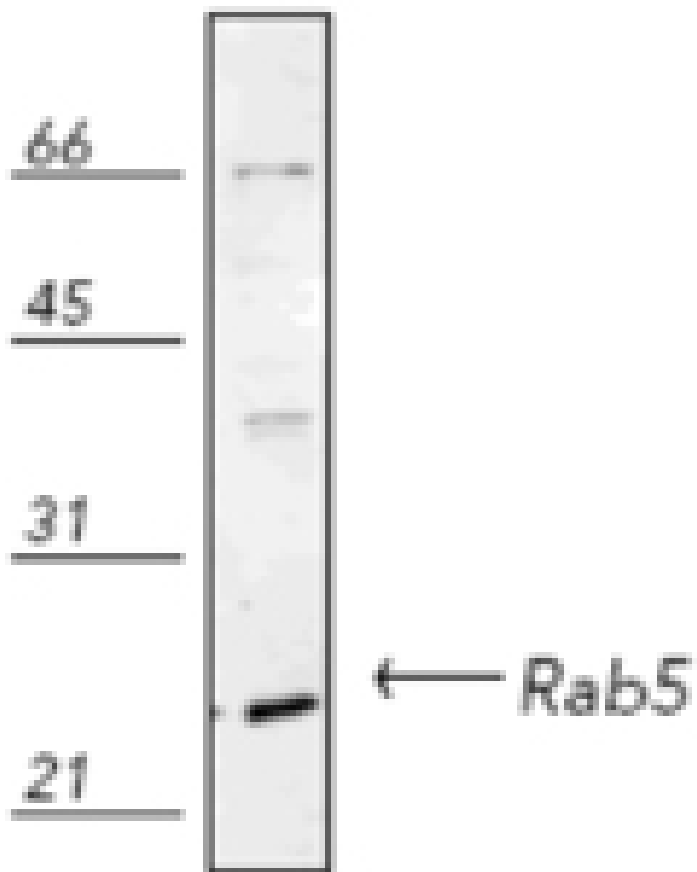
Ordering Information

[Order Online »](#)

ADI-KAP-GP006-E	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of human HeLa (heat shocked) cell lysate, probed with Rab5 pAb.

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 Ras-related protein Rab-5

Application ICC, IHC, WB

Application Notes Detects a band of ~26kDa by Western blot.

Formulation Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.

GenBank ID M35520 (dog)

Host Rabbit

Immunogen Synthetic peptide corresponding to a portion of human Rab5. The sequence is completely conserved in dog.

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Western Blot (1:1,000, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Source Purified from rabbit serum.

Species Reactivity Bovine, Dog, Guinea pig, Hamster, Human, Monkey, Mouse, Rat

UniProt ID P20339 (human), P18066 (dog)

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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)