

RAB4 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: 2

[View Online »](#)

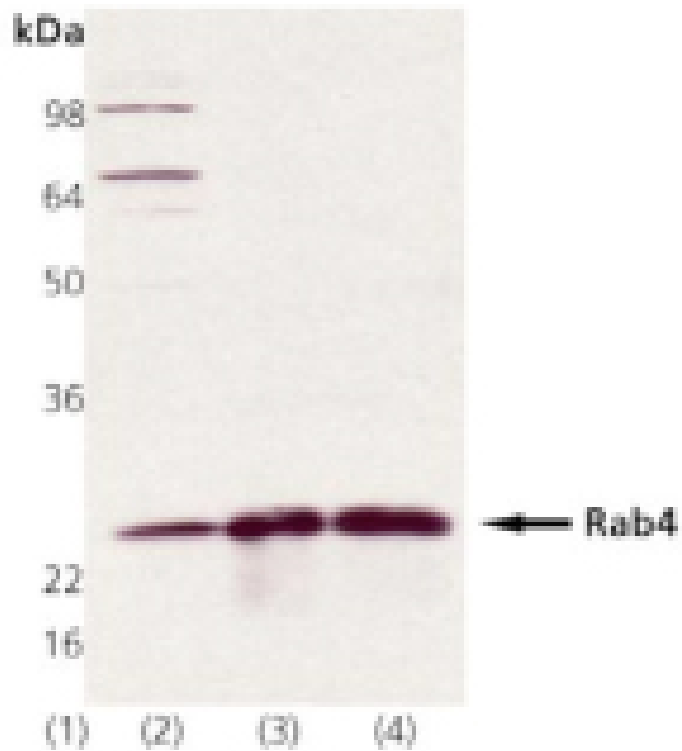
Ordering Information

[Order Online »](#)

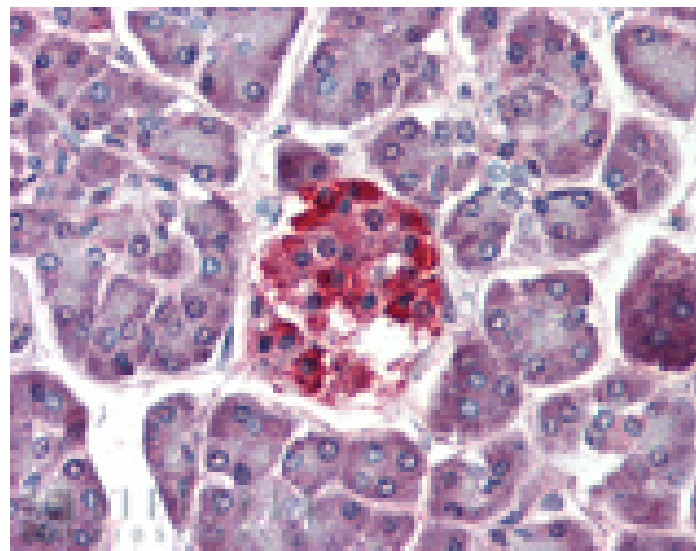
ADI-KAP-GP005-F	200µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Rab4: Lane 1: MW marker, Lane 2: HeLa Cell Lysate (Heat Shocked), Lane 3: Rat Brain Tissue Extract, Lane 4: Mouse Brain Tissue Extract.



Immunohistochemistry analysis of human pancreas tissue stained with RAB4 pAb at 15 μ g/ml.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Ras-related protein Rab-4
Application	ICC, IHC (PS), WB
Application Notes	Detects a band of ~25kDa by Western blot.
Crossreactivity	Weakly cross-reacts with human RAB4.
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.
Host	Rabbit
Immunogen	Recombinant human Rab4.
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:500)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	Mouse, Rat
UniProt ID	P20338

Worry-free Guarantee

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

.



Last modified: May 29, 2024

ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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