

Rap1 polyclonal antibody

The Ras superfamily of small guanosine triphosphatases (GTPases) is traditionally subdivided into five families: Ras, Rho, Rab, Ran, and Arf. Rap1/Krev1 is a member of the Ras family of low molecular weight GTP-binding proteins¹. Ras-like GTPases are ubiquitously expressed, evolutionarily conserved molecular switches that couple extracellular signals to various cellular responses. Rap1 is primarily found at the cytosolic side of intracellular membranes and has two isoforms: Rap1a and 1b2. Both isoforms have a molecular mass of 21 kDa and are isoprenylated at the carboxyl-terminal and phosphorylated by the cAMP-dependent protein kinase A (PKA)³. Rap1 cycles between a GTP-bound active form and a GDP-bound inactive form that is mediated by GTPase activating protein (GAP) and GDP dissociation stimulator (GDS)^{4,5}. Activation occurs by a variety of extracellular stimuli through several conserved guanine nucleotide exchange factors (GEFs) and GTPase activating proteins (GAPs). Rap1 is proposed to regulate Ras mediated signaling and may also be involved in the regulation of integrin-mediated cell adhesion, although the mechanism of regulation is not known. Overexpression of Rap reverses the transformed phenotype induced by Ras, possibly by competing with Ras for interaction with Ras-GAP. Rap has been shown to participate in MAP kinase cascade activated by growth factor and maintaining human T cell anergic state by blocking IL-2 expression.

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

Ordering Information

[Order Online »](#)

ADI-KAP-GP126-D	50µg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Rap1 pAb: Lane 1: Rat Brain Tissue Extract, Lane 2: Mouse Brain Tissue Extract.

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 KREV1

Application WB

Application Notes Detects a band of ~25kDa by Western blot.

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

Host Rabbit

Immunogen Synthetic peptide corresponding to the sequence near the C-terminus of human, mouse and rat Rap1.

Purity Detail Peptide 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, Human, Mouse, Rabbit, Rat

UniProt ID P62834 (Rap1A), P61224 (Rap1B)

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-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

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

Last modified: May 29, 2024