

Rap1 polyclonal antibody

Rap1, which is a member of the Ras family of GTP-binding proteins, cycles between an active GTP-bound and an inactive GDP-bound form that is mediated by GTPase activating protein (GAP). 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.

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

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-KAP-GP125-E	100µg
-----------------	-------

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 KREV1

Application WB

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

Formulation Liquid. In PBS.

GenBank ID A08691

Host Rabbit

Immunogen Synthetic peptide corresponding to a portion of human Rap1a. The sequence is completely conserved in mouse, rat, bovine, chicken, Xenopus. Differs from a portion of Rap1b by 2 amino acids.

Purity Detail Peptide affinity purified.

Recommendation Dilutions/Conditions Western Blot (1µg/ml, 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, Chicken, Dog, Guinea pig, Hamster, Human, Monkey, Mouse, Porcine, Rabbit, Rat, Sheep, Xenopus

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-
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)