

Ras polyclonal antibody

Monomeric G-protein Ras functions as a molecular switch linking receptor and nonreceptor tyrosine kinase activation to a number of downstream signaling pathways. Oncogenic mutations, which occur in approximately 30% of all cases of human cancer, lead to the stimulation of proliferation and the inhibition of apoptosis.

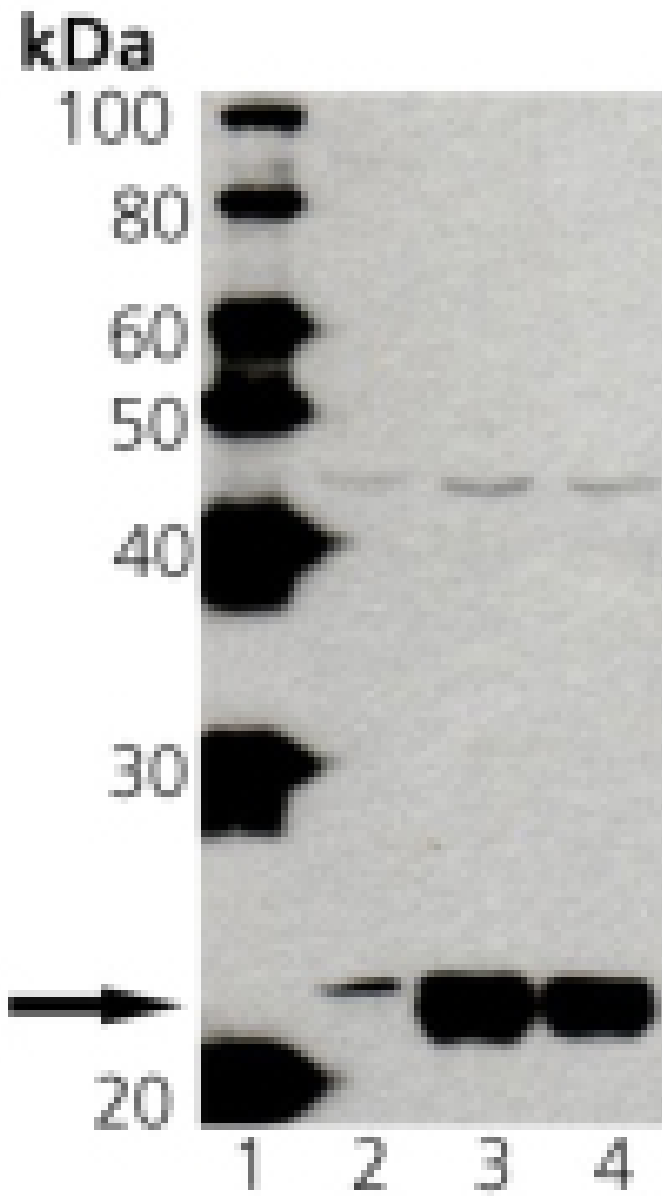
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of anti-Ras pAb: Lane 1: MW marker, Lane 2: HeLa lysate, Mouse Brain Lysate, Rat Brain Lysate.

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application	WB
Application Notes	Detects a band of 21-23kDa by Western blot.
Formulation	Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.
GenBank ID	J00277
Host	Rabbit
Immunogen	Synthetic peptide corresponding to a portion of human Ras.
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:500, 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, Human, Mouse, Rat
UniProt ID	P01112

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

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

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