

# 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 »](#)

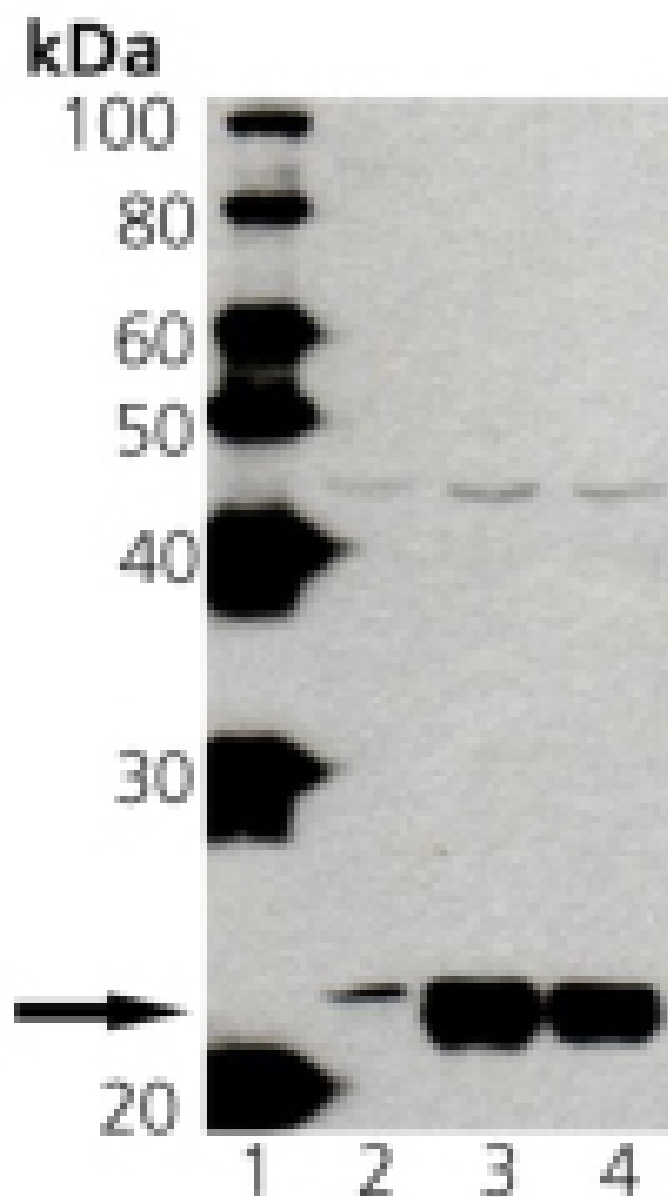
## Ordering Information

[Order Online »](#)

|                 |       |
|-----------------|-------|
| ADI-KAP-GP001-D | 50µg  |
| ADI-KAP-GP001-F | 200µg |

## 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-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)