# 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 »** 

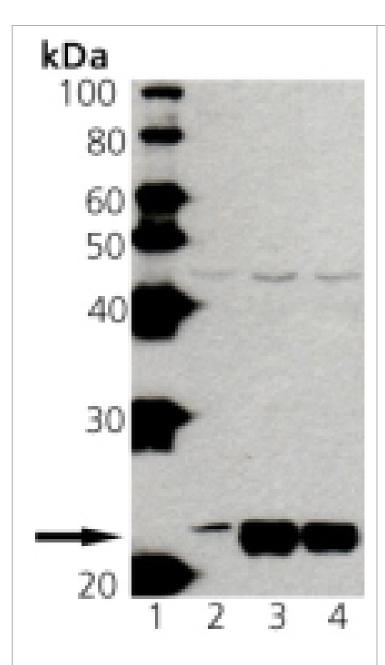
## **Ordering Information**

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



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 m infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

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