
KRAS Polyclonal Antibody, FITC Conjugated

(Catalog #A52270)

Background

Ras proteins bind GDP/GTP and possess intrinsic GTPase activity.

Description

KRAS Polyclonal Antibody, FITC Conjugated. FITC. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human

Isotype

IgG

Uniprot ID

P01116

Purification

>95%, Protein G purified

Immunogen

Recombinant Human GTPase KRas protein (2-168AA)

Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

K-Ras 2, Ki-Ras, c-K-ras, c-Ki-ras, KRAS, KRAS2, RASK2

Application

ELISA; Recommended dilution: ELISA 1:100-1:500
