

RAIDD Polyclonal Antibody (AT100)

(Catalog No. A-0500)

Background

RAIDD, or RIP-associated ICH-1 homologous protein with a death domain, is an adaptor molecule specific for caspase-2 and FASL/TNF receptor-interacting protein RIP. In the presence of RIP and TRADD, CRADD recruits caspase-2 to the TNFR-1 signaling complex.

Description

Rabbit polyclonal antibody to RAIDD

Formulation

Liquid. In PBS containing 0.02% sodium azide.

Specificity

Human

UniProt ID

P78560

Purification

Affinity Purified

Immunogen

Recombinant human RAIDD (aa 1-199)

Storage

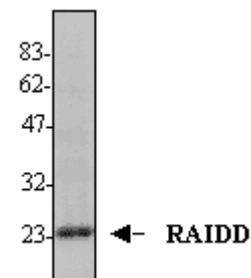
Store at 4°C for short term. Aliquot and store at -20°C for long term. Avoid multiple freeze/thaw cycles.

Alternative Names

CRADD, Caspase and RIP adaptor with death domain, RIP-associated protein with a death domain

Application

IP, WB (1:1000)



▲ Detection of endogenous RAIDD (human) in HEK293T cells by the RAIDD Polyclonal Antibody (AT100).

Ordering Information

Products

RAIDD Polyclonal Antibody (AT100)

Size

50 µg

Cat. No.

A-0500-050

This product is for research purposes only. Not intended for use in diagnostic procedures.