
NAPRT Polyclonal Antibody, HRP Conjugated

(Catalog # A53725)

Background

Catalyzes the conversion of nicotinic acid (NA) to NA mononucleotide (NaMN). Essential for NA to increase cellular NAD levels and prevent oxidative stress of the cells.

Description

NAPRT Polyclonal Antibody, HRP Conjugated. HRP. Raised in: Rabbit.

Formulation

Preservative: 0.1mol/L NaCl, 15mmol/L NaN₃, PH7.2

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q6XQN6

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Nicotinate phosphoribosyltransferase protein (1-310AA)

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

FHA-HIT-interacting protein Nicotinate phosphoribosyltransferase domain-containing protein 1 NAPRT1 FHIP

Application

ELISA, WB; Recommended dilution: ELISA 1:500 - 1:1000, WB 1:1000 – 1:5000
