
GAPDH Polyclonal Antibody, HRP Conjugated

(Catalog #A52539)

Background

Has both glyceraldehyde-3-phosphate dehydrogenase and nitrosylase activities, thereby playing a role in glycolysis and nuclear functions, respectively. Glyceraldehyde-3-phosphate dehydrogenase is a key enzyme in glycolysis that catalyzes the first step of the pathway by converting D-glyceraldehyde 3-phosphate (G3P) into 3-phospho-D-glyceroyl phosphate. Modulates the organization and assembly of the cytoskeleton. Also participates in nuclear events including transcription, RNA transport, DNA replication and apoptosis. Nuclear functions are probably due to the nitrosylase activity that mediates cysteine S-nitrosylation of nuclear target proteins

Description

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

Formulation

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

Specificity

Chicken

Isotype

IgG

Uniprot ID

P00356

Purification

Protein G purified

Immunogen

Recombinant Chicken Glyceraldehyde-3-phosphate dehydrogenase protein (2-333AA)

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

Peptidyl-cysteine S-nitrosylase GAPDH, GAPDH, GAPD

Application

ELISA; Recommended dilution: ELISA 1:500-1:1000
