
GAPDH Tag Antibody

(Catalog No. A26005)

Description

GAPDH Tag Antibody recognizes glyceraldehyde-3-phosphate dehydrogenase (GAPDH), a 36kD multifunctional protein whose main function is to catalyse the reversible oxidative phosphorylation of glyceraldehyde-3-phosphate, in conjunction with inorganic phosphate and nicotinamide adenine dinucleotide (NAD). This reaction is an important energy yielding step in carbohydrate metabolism. GAPDH has also been shown to translocate to the nucleus under a variety of stressors, most of which are associated with oxidative stress, whereby it mediates cell death. A further report has shown that GAPDH binds to several proteins that are responsible for neurodegenerative diseases, such as amyloid precursor protein and Huntingtin.

Antibody Type

Monoclonal Antibody

Purification

Purified IgG - liquid

Immunogen

Human cardiac muscle GAPDH

Isotype

IgG1

Size

200 µg

Formulation

Phosphate buffered saline. 0.09% Sodium Azide (NaN₃)

Specificity

Bovine; Cat; Dog; Fish; Goat; Mouse; Pig; Rabbit; Rat

Storage

Store at +4°C or at -20°C if preferred. Storage in frost-free freezers is not recommended. This product should be stored undiluted. Avoid repeated freezing and thawing as this may denature the antibody.

Handling Recommendations

Should this product contain a precipitate we recommend microcentrifugation before use. For maximum recovery of the products, centrifuge the vial prior to opening the cap.

Applications & Suggested Dilutions

ELISA; Immunofluorescence; Immunohistology - Frozen; Immunoprecipitation; Western Blotting

Ordering Information:

Products	Size	Cat. No.
GAPDH Tag Antibody	200 µg	A26005-1

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