

NME5 Polyclonal Antibody

(Catalog # A63030)

Background

Does not seem to have NDK kinase activity. Confers protection from cell death by Bax and alters the cellular levels of several antioxidant enzymes including Gpx5. May play a role in spermiogenesis by increasing the ability of late-stage spermatids to eliminate reactive oxygen species.

Description

NME5 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

P56597

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Nucleoside diphosphate kinase homolog 5 protein (1-212AA)

Storage

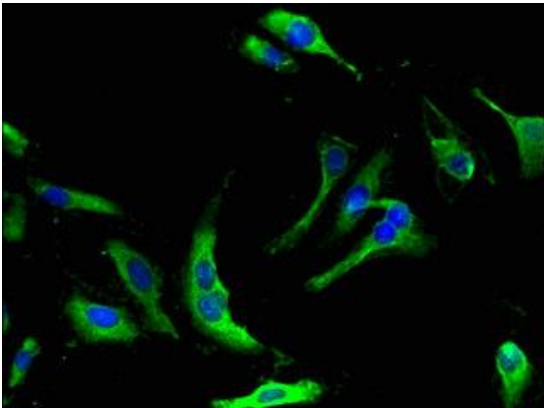
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Nucleoside diphosphate kinase homolog 5, NDK-H 5, NDP kinase homolog 5, Inhibitor of p53-induced apoptosis-beta, IPIA-beta, Testis-specific nm23 homolog, nm23-H5, NME5

Application

ELISA, IF; Recommended dilution: IF:1:50-1:200



Immunofluorescent analysis of HeLa cells at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)