

PAX3 Polyclonal Antibody

(Catalog # A50404)

Background

Transcription factor that may regulate cell proliferation, migration and apoptosis. Involved in neural development and myogenesis.

Description

PAX3 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

P23760

Purification

Protein G purified

Immunogen

Recombinant Human Paired box protein Pax-3 protein (90-184AA)

Storage

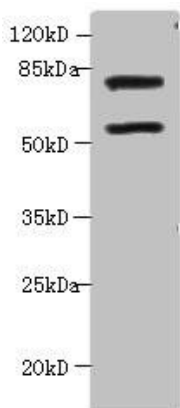
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Paired box protein Pax-3, HuP2, PAX3, HUP2

Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:5000, IHC:1:20-1:200, IF:1:50-1:200



Western blot

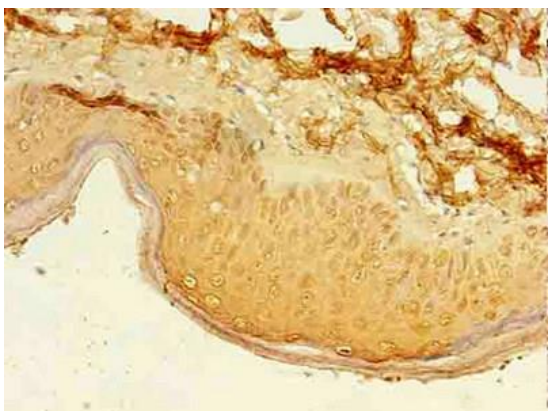
All lanes: PAX3 Polyclonal Antibody at 2ug/ml+Hela cells

Secondary

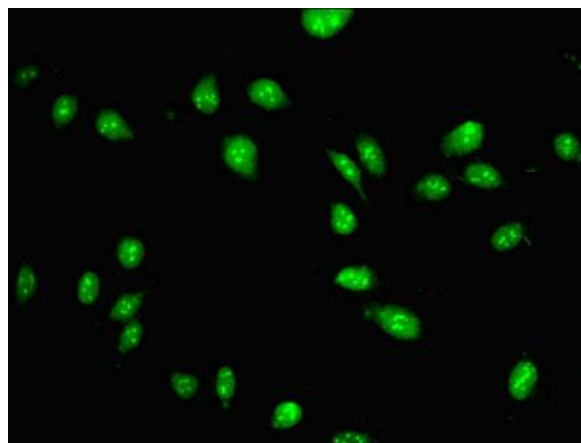
Goat polyclonal to rabbit at 1/10000 dilution

Predicted band size: 53, 25, 23, 45, 46, 54, 54, 56 kDa

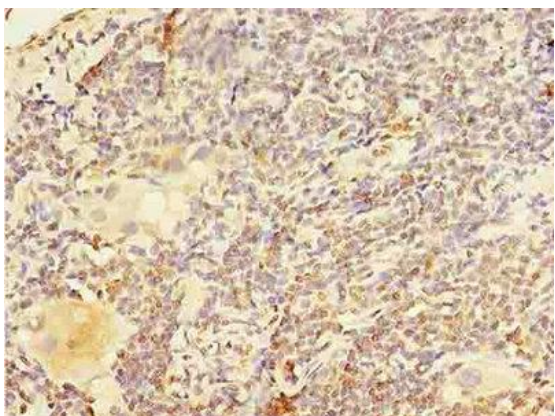
Observed band size: 53 kDa



Immunohistochemistry of paraffin-embedded human skin using PAX3 Polyclonal Antibody at dilution 1:100



Immunofluorescent analysis of Hela cells using PAX3 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)



Immunohistochemistry of paraffin-embedded human lung cancer using PAX3 Polyclonal Antibody at dilution 1:100