

PADI3 Polyclonal Antibody

(Catalog # A68054)

Background

Catalyzes the deimination of arginine residues of proteins.

Description

PADI3 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

Q9ULW8

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Protein-arginine deiminase type-3 protein (1-664AA)

Storage

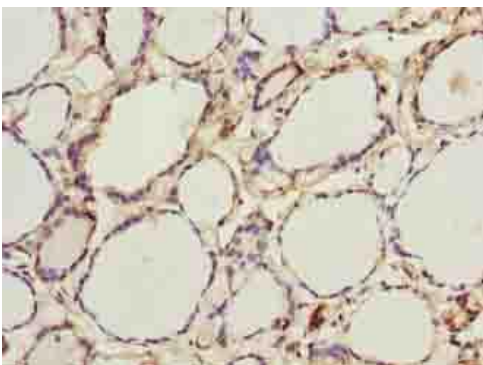
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

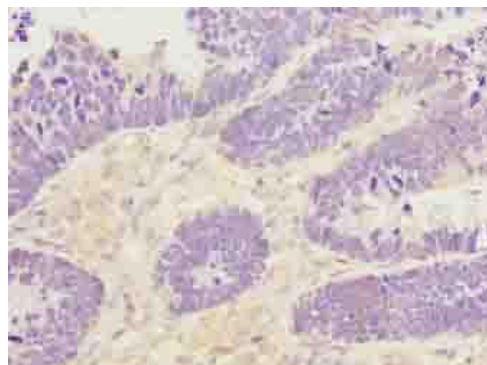
Protein-arginine deiminase type-3, Peptidylarginine deiminase III, Protein-arginine deiminase type III, PADI3, PAD3, PDI3

Application

ELISA, IHC; Recommended dilution: IHC:1:20-1:200



Immunohistochemistry of paraffin-embedded human thyroid tissue using PADI3 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human ovarian cancer using PADI3 Polyclonal Antibody at dilution of 1:100