
NID1 Polyclonal Antibody, FITC Conjugated

(Catalog #A68637)

Background

Sulfated glycoprotein widely distributed in basement membranes and tightly associated with laminin. Also binds to collagen IV and perlecan. It probably has a role in cell-extracellular matrix interactions.

Description

NID1 Polyclonal Antibody, FITC Conjugated. FITC. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

P14543

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Nidogen-1 protein (792-942AA)

Storage

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

Alternative Names

Nidogen-1, NID-1, Entactin, NID1, NID

Application

ELISA; Recommended dilution: ELISA 1:100-1:500
