
LAMA5 Polyclonal Antibody

(Catalog #A59602)

Background

Binding to cells via a high affinity receptor, laminin is thought to mediate the attachment, migration and organization of cells into tissues during embryonic development by interacting with other extracellular matrix components.

Description

LAMA5 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

O15230

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Laminin subunit alpha-5 protein (3401-3692AA)

Storage

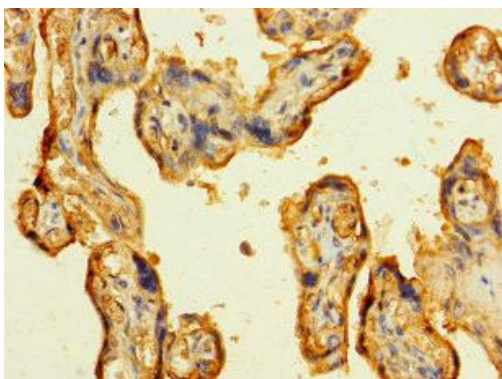
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Laminin subunit alpha-5, LAMA5, KIAA0533, KIAA1907, Laminin-10 subunit alpha, Laminin-11 subunit alpha, Laminin-15 subunit alpha

Application

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



Immunohistochemistry of paraffin-embedded human placenta tissue using LAMA5 Antibody at dilution of 1:100