
KLK5 Polyclonal Antibody

(Catalog #A51429)

Background

May be involved in desquamation.

Description

KLK5 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

Q9Y337

Purification

Protein G purified

Immunogen

Recombinant Human Kallikrein-5 protein (23-293AA)

Storage

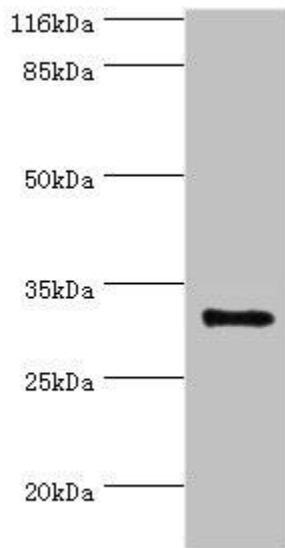
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

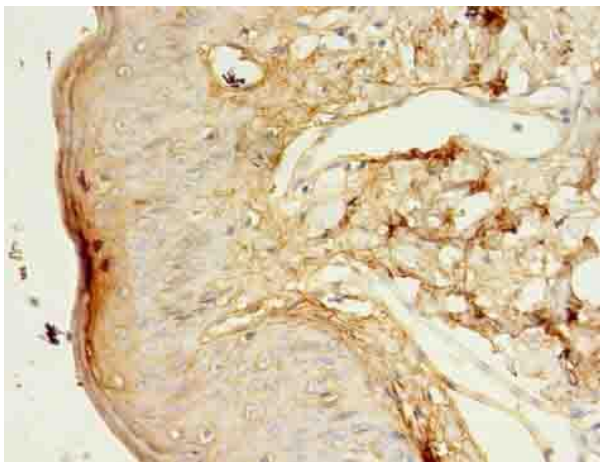
Kallikrein-5, Kallikrein-like protein 2, KLK-L2, Stratum corneum tryptic enzyme, KLK5, SCTE

Application

ELISA, WB, IHC; Recommended dilution: WB:1:500-1:2000, IHC:1:20-1:200



Western blot
All lanes: KLK5 Polyclonal Antibody at 2ug/ml+MCF-7
whole cell lysate
Secondary
Goat polyclonal to rabbit at 1/10000 dilution
Predicted band size: 32kDa
Observed band size: 32kDa



Immunohistochemistry of paraffin-embedded human skin
tissue using KLK5 Polyclonal Antibody at dilution of
1:100