
Klk1b22 Polyclonal Antibody

(Catalog A59578)

Background

Glandular kallikreins cleave Met-Lys and Arg-Ser bonds in kininogen to release Lys-bradykinin.

Description

Klk1b22 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Mouse

Isotype

IgG

Uniprot ID

P15948

Purification

>95%, Protein G purified

Immunogen

Recombinant Mouse Kallikrein 1-related peptidase b22 protein (25-259AA)

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 1-related peptidase b22 (EC:3.4.21.35), Klk1b22, Klk-22, Klk22, Beta-NGF-endopeptidase, Epidermal growth factor-binding protein type A, EGF-BP A, Glandular kallikrein K22, mGK-22, Nerve growth factor beta chain endopeptidase, Tissue kallikrein 22

Application

ELISA; Recommended dilution: ELISA 1:500-1:2000
