
PLAT Polyclonal Antibody, HRP Conjugated

(Catalog #A56468)

Background

Converts the abundant, but inactive, zymogen plasminogen to plasmin by hydrolyzing a single Arg-Val bond in plasminogen. By controlling plasmin-mediated proteolysis, it plays an important role in tissue remodeling and degradation, in cell migration and many other physiopathological events. Plays a direct role in facilitating neuronal migration.

Description

PLAT Polyclonal Antibody, HRP Conjugated. HRP. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

P00750

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Tissue-type plasminogen activator protein (36-556AA)

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

t-plasminogen activator, PLAT

Application

ELISA; Recommended dilution: ELISA 1:500-1:1000
