

Thpo Polyclonal Antibody

(Catalog # A64108)

Background

Lineage-specific cytokine affecting the proliferation and maturation of megakaryocytes from their committed progenitor cells. It acts at a late stage of megakaryocyte development. It may be the major physiological regulator of circulating platelets.

Description

Thpo Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Rat, Mouse, Human

Isotype

IgG

Uniprot ID

P49745

Purification

Protein G purified

Immunogen

Recombinant Rat Thrombopoietin protein (22-194AA)

Storage

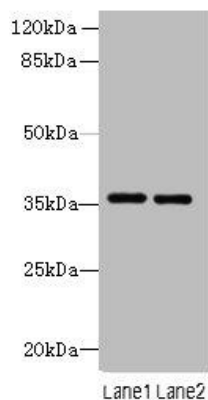
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Thrombopoietin, Thpo

Application

ELISA, WB; Recommended dilution: WB:1:1000-1:5000



Western blot

All lanes: Thpo Polyclonal Antibody at 4ug/ml

Lane 1: Mouse liver tissue

Lane 2: 293T whole cell lysate

Secondary

Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 35 kDa

Observed band size: 35 kDa