

igg Light Chain Polyclonal Antibody

(Catalog # A57793)

Description

igg Light Chain Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Purification

>95%, Protein G purified

Immunogen

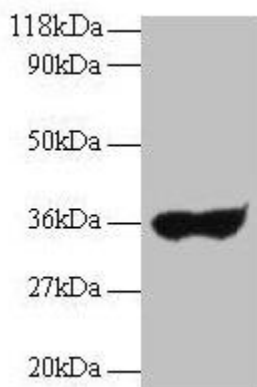
Human IgG light chain protein (Native Protein)

Storage

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

Application

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



Western blot

All lanes: Human Igg light chain antibody at 2ug/ml + Human serum at 1: 100

Secondary

Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 35 kDa

Observed band size: 35 kDa