
EPO Polyclonal Antibody, HRP Conjugated

(Catalog # A64093)

Description

EPO 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

P01588

Purification

Protein G purified

Immunogen

Recombinant Human Erythropoietin protein (28-191AA)

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

Erythropoietin, Epoetin, EPO

Application

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