

PYC1 Polyclonal Antibody

(Catalog # A53675)

Background

Pyruvate carboxylase catalyzes a 2-step reaction, involving the ATP-dependent carboxylation of the covalently attached biotin in the first step and the transfer of the carboxyl group to pyruvate in the second.

Description

PYC1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Baker's yeast

Isotype

IgG

Uniprot ID

P11154

Purification

>95%, Protein G purified

Immunogen

Recombinant Baker's yeast Pyruvate carboxylase 1 protein (18-470AA)

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

PYC1, PYV, YGL062W

Application

ELISA
