

Cytochrome P450 Polyclonal Antibody

(Catalog No. A-0482)

Background

Cytochrome p450 is a highly conserved and diverse family of heme-containing proteins that function in the metabolism of numerous endogenous and exogenous substrates. The human genome encodes 57 p450 genes divided among 18 families, which together account for the diversity of substrates metabolized. p450 is the primary conjugating enzyme in phase I metabolism and detoxification of endogenous and exogenous compounds, and are primarily localized to the inner membranes of mitochondria and the ER. Eukaryotic microsomal cytochrome p450 enzymes receive electrons from the FAD- and FMN-containing enzyme NADPH:cytochrome p450 reductase (CPR).

Description

Rabbit polyclonal antibody to Cytochrome P450

Formulation

Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.

Immunogen

Native rat Cytochrome p450.

Specificity

Human, Mouse, Rat, Clam, Dog, Guinea Pig, Horse, Monkey, Rabbit

Purification

Protein A-affinity purified.

Storage

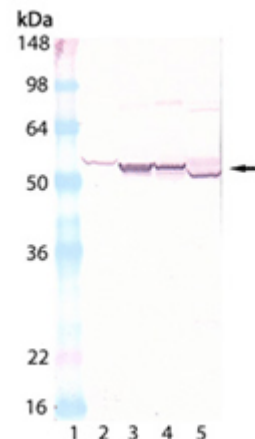
Store at -20°C. Avoid multiple freeze/thaw cycles.

Alternative Names

CYPB

Application

WB (1:1000)



▲Western blot analysis of Cytochrome P450: Lane 1: MWM, Lane 2: Human Liver microsomes, Lane 3: Rat Liver Microsomes, Lane 4: Mouse Liver Microsomes, Lane 5: Rabbit Liver Microsomes.

Ordering Information

Products

Cytochrome P450 Polyclonal Antibody

Size

200 µg

Cat. No.

A-0482-200

This product is for research purposes only. Not intended for use in diagnostic procedures.