Cytochrome P450 3A1 (rat) polyclonal antibody

Members of the cytochrome P-450 3A (CYP3A) sub-family catalyze the 6β-hydroxylation of testosterone, and the metabolism of a number of drugs such as cyclosporin A (Prod. No. BML-A195), and warfarin. They are inducible by glucocorticoids such as dexamethasome (Prod. No. BML-EI126) and by macrolide antibiotics such as erythromycin and pregnenolone 16α-carbonitrile (U-14975, PCN) (Prod. No. BML-S620). Rat CYP3A1 and CPY3A2 share 89% sequence similarity at the amino acid level and are predominantly located in the liver, principally in the centrilobular region. PCN induces CYP3A1 gene expresion in regions of the intestine, lung and kidney.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

BML-CR3315-0025 25µl

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name CYP3A1

Application IHC (FS), IHC (PS), WB

Formulation Liquid. Serum containing 0.01% sodium azide.

Host Sheep

Immunogen Synthetic peptide corresponding to a portion of C-terminal

rat CYP3A1.

Species Reactivity Rat

Specificity Recognizes rat CYP3A1. Not suitable for use on human

tissue.

UniProt ID P04800

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

Last modified: May 29, 2024

