

Cytochrome P450 3A1 (rat) polyclonal antibody

Members of the cytochrome P-450 3A (CYP3A) sub-family catalyze the 6 β -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 dexamethasone (Prod. No. BML-EI126) and by macrolide antibiotics such as erythromycin and pregnenolone 16 α -carbonitrile (U-14975, PCN) (Prod. No. BML-S620). Rat CYP3A1 and CYP3A2 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 expression in regions of the intestine, lung and kidney.

This antibody is covered by our [Worry-Free Guarantee](#).

Ordering Information

[Order Online »](#)

BML-CR3315-0025	25 μ l
-----------------	------------

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](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)