

Cytochrome P450 2D4 (rat) polyclonal antibody

Cytochrome P-450 2D4 (CYP2D4; also known as debrisoquine-4-hydroxylase), is found at very low levels in normal rat brain, but administration of clozapine (Prod. No. BML-D107) induces high levels in many areas including the *substantia nigra pars compacta*, ventral tegmental area, in the olfactory bulb and in granule and Purkinje cells in the cerebellum.

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

Ordering Information

[Order Online »](#)

BML-CR3240-0025	25µl
BML-CR3240-0100	100µl

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 CYP2D4

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

Formulation Liquid. Serum containing 0.01% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to the C-terminus of rat liver CYP2D4.

Species Reactivity Rat

Specificity Recognizes rat CYP2D4. Not suitable for use on human tissue.

UniProt ID P13108

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)