

# 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 <a href="#">Worry-Free Guarantee</a>

Last modified: May 29, 2024



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)