

# Cytochrome P450 1A1 (rat) polyclonal antibody

Members of the cytochrome P-450 1A (CYP1A) sub-family are induced by a range of polycyclic aromatic compounds such as benzpyrene, 3-methylcholanthrene (Prod. No. BML-GR239), and *pro*-carcinogenic heterocyclic aromatic amines. Such compounds have been reported to induce Cytochrome P-450 (CYP1A1) in rat liver, human placenta and lung. A rare polymorphism of the CYP1A1 gene has been implicated in familial Parkinson's Disease (PD). In the rat, CYP1A1 has been localized to the *globus pallidus*, but not to the *substantia nigra* where cell death associated with the pathology of PD occurs.

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

## Ordering Information

[Order Online »](#)

BML-CR3110-0025	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 CYP1A1

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

Formulation Liquid. Serum containing 0.01% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to a portion of C-terminal rat CYP1A1.

Species Reactivity Rat

Specificity Recognizes rat CYP1A1. Not suitable for use on human tissue.

UniProt ID P00185

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)