

Cytochrome P450 1A2 (human) 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. Cigarette smoking has been reported to induce CYP1A2 in human liver, and appears to be a major catalyst in the hepatic activation of PhIP, (2-amino-1-methyl-6-phenylimidazo[4,5,b]-pyridine). The mono-oxygenase inhibitor enoxacin induces CYP1A2 in rat liver, whereas a related compound ciprofloxacin does not.

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

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-CR3130-0025

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	CYP1A2
Application	IHC (FS), IHC (PS), WB
Formulation	Liquid. Serum containing 0.01% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to the internal domain (aa 284-298) of human CYP1A2.
Species Reactivity	Human
Specificity	Recognizes human CYP1A1. Not suitable for use on rat tissue.
UniProt ID	P05177
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee

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-
fr@enzolifesciences.com](mailto:info-fr@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)