

# Cytochrome P450

## polyclonal antibody

Cytochrome p450 is a highly conserved and diverse family of heme-containing proteins that function in the metabolism of numerous endogenous and exogenous substrates.. The human genome encodes 57 p450 genes divided among 18 families, which together account for the diversity of substrates metabolized. p450 is the primary conjugating enzyme in phase I metabolism and detoxification of endogenous and exogenous compounds, and are primarily localized to the inner membranes of mitochondria and the ER. Eukaryotic microsomal cytochrome p450 enzymes receive electrons from the FAD- and FMN-containing enzyme NADPH:cytochrome p450 reductase (CPR).

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

Citations: 5

[View Online »](#)

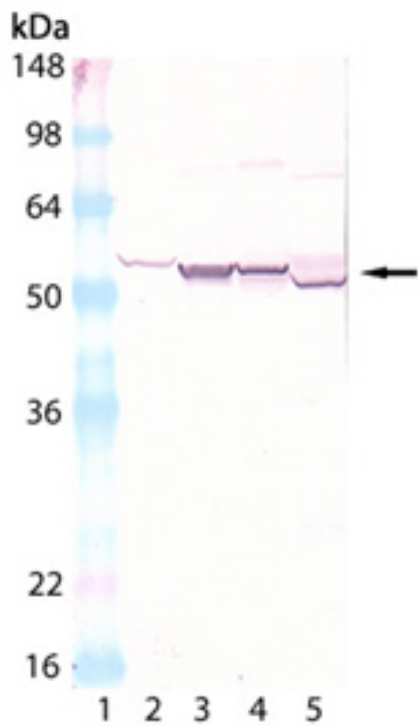
### Ordering Information

[Order Online »](#)

ADI-MFO-100-D	50µg
ADI-MFO-100-F	200µg

### Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Cytochrome P450: Lane 1: MWM, Lane 2: Human Liver microsomes, Lane 3: Rat Liver Microsomes, Lane 4: Mouse Liver Microsomes, Lane 5: Rabbit Liver Microsomes.

<b>Handling</b>	Avoid freeze/thaw cycles.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

## Product Details

Alternative Name	CYPR
Application	WB
Application Notes	Detects a band of ~50-55 kDa by Western blot.
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.
GenBank ID	M20131
Host	Rabbit
Immunogen	Native rat Cytochrome p450.
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from rabbit serum.
Species Reactivity	Clam, Dog, Guinea pig, Horse, Human, Monkey, Mouse, Rabbit, Rat
UniProt ID	P05182
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>



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)