

Cytochrome P450 2C8, 2C9, 2C19 (human) and 2C12 (rat) polyclonal antibody

Mephenytoin 4-hydroxylase family of CYPs. Metabolizes anti-depressants.
Cytochrome P-450 (CYP2C19) is a sex-dependent steroid hydroxylase.

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

Ordering Information

[Order Online »](#)

BML-CR3280-0025	25µl
BML-CR3280-0100	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 CYP2C8, 2C9, 2C19 and CYP2C12

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 human liver CYP2C8, 2C9, 2C19, and rat 2C12.

Species Reactivity Human

Specificity Recognizes human CYP2C8, CYP2C9, CYP2C19 and rat CYP2C12.

UniProt ID P10632 (human CYP2C8), P11712 (human CYP2C9), P33261 (human CYP2C19), P11510 (rat CYP2C12)

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)