

ACE1 (human) (CT) polyclonal antibody

ACE1 (angiotensin converting enzyme-1, ACE) is a key regulator of blood pressure.

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

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	Angiotensin converting enzyme 1, ACE
Application	WB
Crossreactivity	Does not cross-react with ACE2.
Formulation	Liquid. In PBS containing 0.05% sodium azide and 50% glycerol.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to the second catalytic domain (C-domain) of the human ACE1.
Species Reactivity	Human
Specificity	Recognizes somatic and testicular ACE1 and their C-terminal fragments.
UniProt ID	P12821
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-
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)