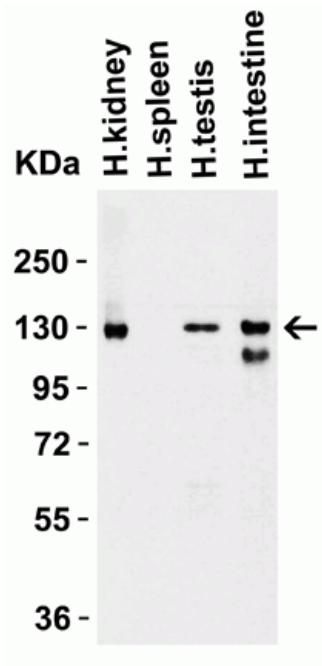


ACE2 polyclonal antibody

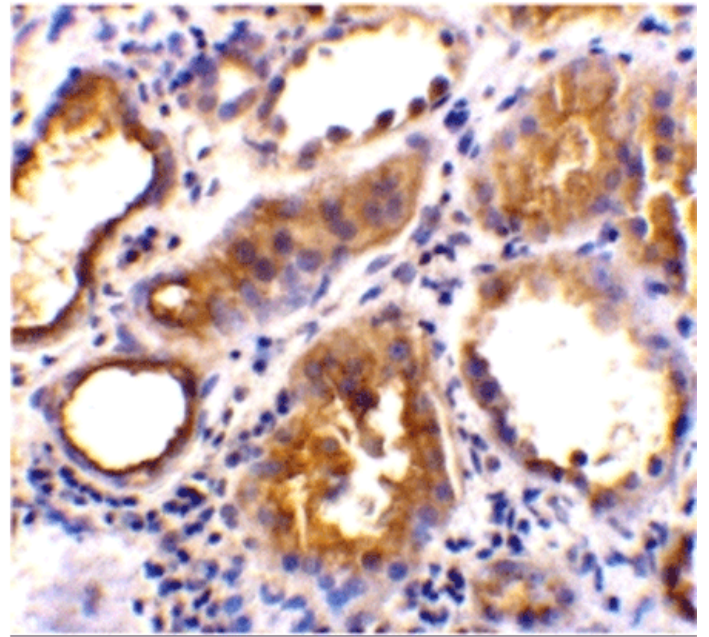
This antibody is covered by our [Worry-Free Guarantee](#).

Manuals, SDS & CofA

[View Online »](#)



Western Blot validation in Human Tissues



Immunohistochemical analysis of paraffin-embedded human kidney tissue using ACE2 antibody at 2 $\mu\text{g/ml}$.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Angiotensin converting enzyme 2, SARS-CoV receptor, SARS-CoV2 receptor

Application ELISA, IF, IHC, WB

Formulation Liquid. In PBS containing 0.02% sodium azide.

Host Rabbit

Immunogen Peptide corresponding to 17 amino acids near the center of human ACE2. The immunogen is located within amino acids 180 – 230 of ACE2.

Isotype IgG

Purity Detail Affinity purified.

Species Reactivity Human, Mouse, Rat

UniProt ID Q9BYF1

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)