

Polyclonal Antibody to ACE2 (IN1)

Ordering Information

[Order Online »](#)

PSC-3215-C100	100µg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	1 year.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Angiotensin-converting Enzyme 2, SARS Receptor
Application	WB
Crossreactivity	No cross response to ACE1.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 192-208 of human ACE2 (Angiotensin-converting enzyme 2).
Positive Control	Prod. No. PSC-1305, Kidney Cell Lysate
Species Reactivity	Human, Mouse
Technical Info / Product Notes	For the Original Manufacturer's data sheet please click here .



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com