

Angiotensin II Converting Enzyme Activity Assay

A plate-based fluorometric assay to measure Angiotensin II converting enzyme (ACE2) in

various samples.

Angiotensin II converting enzyme (ACE2) is part of the renin-angiotensin system (RAS). It plays a pivotal role in both blood pressure regulation and entry of coronavirus. It cleaves Angiotensin II converting it into Angiotensin 1-7. Whereas Angiotensin II is a vasoconstrictor, Angiotensin 1-7 is a vasodilator, so the action of ACE-2 can exert significant control over blood pressure. ACE2 is a known receptor of human coronaviruses, such as SARS. It is a cell-surface protein, expressed in vascular endothelial cells of lung, kidney and heart. ACE2 is a possible therapeutic target for cardiovascular and coronavirus-induced diseases. AkrivisBio's ACE2 Activity Assay is a simple sensitive means of monitoring ACE-2 activity. We utilize a synthetic peptide substrate which releases MCA upon ACE2 based cleavage. A specific ACE-2 inhibitor is included allowing for differentiation of ACE2 activity from other proteolytic activities which may recognize the peptide substrate.

Ordering Information

[Order Online »](#)

AKR-MA-0158	100 wells
-------------	-----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Fluorescent detection

Species Reactivity Species independent

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