

ANK1 antibody – C-terminal region

This is a rabbit polyclonal antibody against ANK1. It was validated on Western Blot and immunohistochemistry by Aviva Systems Biology.

Ordering Information

[Order Online »](#)

ASB-ARP42566-T100	100µl
-------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application IHC, WB

Host Rabbit

Immunogen The immunogen for anti-ANK1 antibody: synthetic 67 directed towards the C terminal of human ANK1

Reconstitution Add 100 µl of distilled water. Final anti-ANK1 antibody concentration is 1 mg/ml in PBS buffer.

Species Reactivity Human, Mouse, Rat, Xenopus

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).

UniProt ID Q4KMP7



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