

IRF6 Antibody – middle region

Ordering Information

[Order Online »](#)

ASB-ARP31995-P050	100µl
-------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Host Rabbit

Immunogen The immunogen is a synthetic peptide directed towards the middle region of human IRF6

Purity Detail Affinity Purified

Species Reactivity Human

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

UniProt ID O14896



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