

hopQ H. Pylori Antibody

Ordering Information

[Order Online »](#)

FGX-HOPQ-101AP	100µg
----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Regulatory Status RUO - Research Use Only

Product Details

Application	ELISA, ICC, IF, IHC, IP, WB
Host	Rabbit
Immunogen	Heat killed and denatured Helicobacter pylori extracts
Isotype	IgG
Purity Detail	Affinity Purified
Species Reactivity	Helicobacter pylori
UniProt ID	J0TJV0



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