

MURF1 Antibody

Ordering Information

[Order Online »](#)

FGX-MURF1-101AP	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Regulatory Status RUO - Research Use Only

Product Details

Application	ELISA, IHC, IP, WB
Host	Rabbit
Immunogen	Synthetic peptide taken within amino acid region 270-310 on human MURF1 protein.
Isotype	IgG
Purity Detail	Affinity Purified
Species Reactivity	Human, Mouse, Rat
UniProt ID	Q969Q1



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