

Phospho-MERTK Antibody

Ordering Information

[Order Online »](#)

FGX-PMKT-140AP	100µg
----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Regulatory Status RUO - Research Use Only

Product Details

Application	ELISA, Flow Cytometry, ICC, IF, IHC, WB
Host	Rabbit
Immunogen	Synthetic peptide corresponding to amino acids 746-757 of Human MERTK. Sequence:CKKIYpSGDYpYpRQG
Species Reactivity	Human, Monkey, Mouse, Rat
UniProt ID	P57097



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