

TRAIL (human) monoclonal antibody (III6F)

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-326-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored as recommended.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Apo-2L, TNFSF 10, CD253
Application	ELISA, Flow Cytometry, IHC (FS), IP, WB
Clone	III6F
Formulation	Liquid. In PBS containing 0.02% sodium azide.
GenBank ID	U57059
Gene/Protein Identifier	8743 (Entrez GeneID)
Host	Mouse
Immunogen	Recombinant human soluble TRAIL.
Isotype	IgG2b
Source	Purified from hybridoma tissue culture supernatant.
Species Reactivity	Human
UniProt ID	P50591



ENZO LIFE SCIENCES,
INC.
Phone: +1 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(EU) LTD
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
ENZO LIFE SCIENCES
Phone: +32 23 466 0426
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

This antibody is covered by our **Worry-Free Guarantee**