

TRAIL-R2 (human):Fc (human), (recombinant) (FITC conjugate)

Citations: 10

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-522-005F-C050	FITC - 50µg
-------------------	-------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored at -20°C.
Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	DR5, TNFRSF 10B, KILLER, CD262
Application Notes	Flow Cytometry: detection of TRAIL cell surface expression.
Biological Activity	Inhibits human soluble TRAIL-mediated apoptosis.
Concentration	1mg/ml.
Endotoxin Content	<0.1EU/µg purified protein (LAL test; Bio Whittaker).
Formulation	Liquid. In PBS containing 0.02% sodium azide.
MW	~46kDa (SDS-PAGE).
Purity	≥95% (SDS-PAGE)
Source	Produced in HEK 293 cells. The extracellular domain of human TRAIL-R2 (DR5) (aa 52-212) is fused to the Fc portion of human IgG1.
Specificity	Binds human and mouse TRAIL.
Technical Info / Product Notes	Historical data has shown inhibition of human soluble TRAIL-induced apoptosis in a concentration range of 0.5-10µg/ml.
UniProt ID	O14763



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

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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