

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

Citations: 15

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

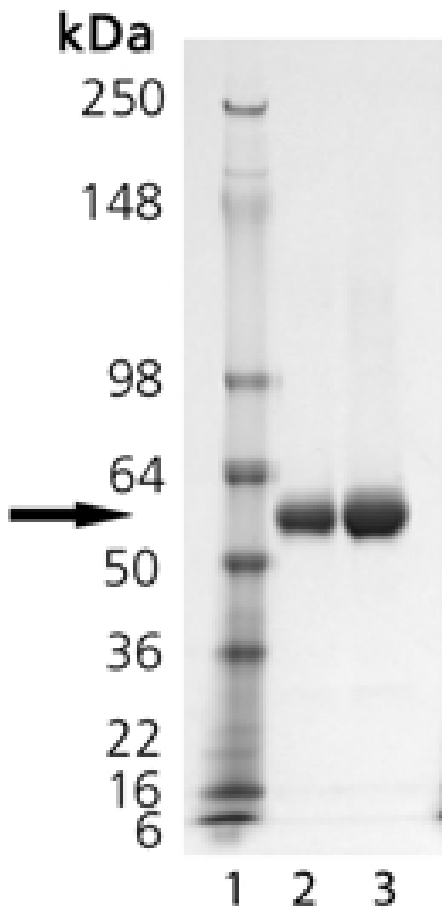


Fig 1: SDS-PAGE Analysis. Lane 1: MW Marker, Lane 2: 1 μ g, Lane 3: 2 μ g TRAIL-R2.

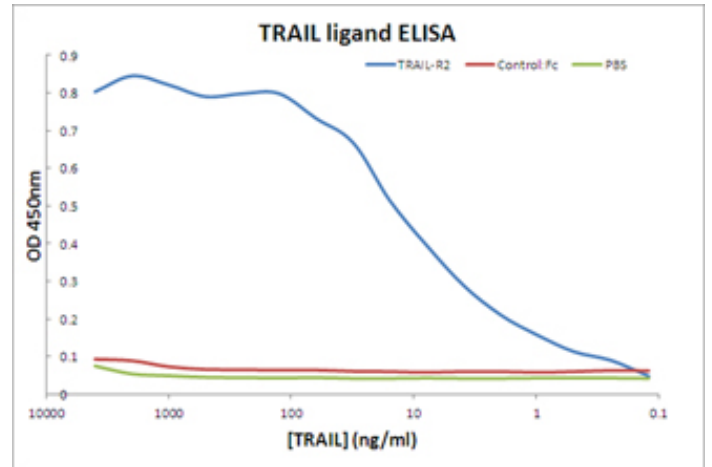


Fig 2: ELISA Analysis: 100 ng/well of receptor or control is coated on a 96-well plate. After blocking, ligand is coated on a 96-well plate. After blocking, ligand is titrated 2-fold in series, and detected with anti-flag antibody. TRAIL (Prod # ALX-522-003) shows positive for binding to TRAIL-R2 and negative for binding to either Fc control protein (Prod # ALX-203-004) and buffer control.

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	ELISA: binds to TRAIL at 10-100 ng/ml.
Biological Activity	Inhibits human soluble TRAIL-mediated apoptosis.
Concentration	1mg/ml after reconstitution.
Endotoxin Content	<0.1EU/μg purified protein (LAL test; Bio Whittaker).
Formulation	Lyophilized. Contains PBS.
MW	58kDa (SDS-PAGE).
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute with 50μl sterile water. Further dilutions should be made with medium containing 5% fetal calf serum or a carrier protein.
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.

Last modified: May 29, 2024



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)