

TRAIL-R1 (human) monoclonal antibody (HS101) (ATTO 647N conjugate)

TRAIL-R1 is a receptor for the cytotoxic ligand TRAIL. The adapter molecule FADD recruits caspase-8 to the activated receptor. The resulting death-inducing signaling complex (DISC) performs caspase-8 proteolytic activation which initiates the subsequent cascade of caspases (aspartate-specific cysteine proteases) mediating apoptosis. Promotes the activation of NF- κ B.

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

Citations: 11

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-297TS-T100

ATTO 647N - 100 tests

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Do not freeze. Protect from light.
Short Term Storage	+4°C
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	TRAIL receptor 1, Apo-2, CD261, DR4, Death receptor 4, TNFRSF 10A, TNF-related apoptosis-inducing ligand receptor 1, Tumor necrosis factor receptor superfamily member 10A
Application	Flow Cytometry, ICC
Clone	HS101
Crossreactivity	Does not cross-react with human TRAIL-R2, -R3 or -R4.
Formulation	Liquid. 100µg in PBS containing 0.02% sodium azide.
GenBank ID	U90875
Gene/Protein Identifier	8797 (Entrez GeneID)
Host	Mouse
Immunogen	Recombinant human TRAIL-R1.
Isotype	IgG1
Recommendation Dilutions/Conditions	Flow Cytometry (1:200)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human

Technical Info / Product Notes

ATTO-fluorescent antibodies show increased photostability, outstanding brightness and intense signals. ATTO dyes are thermally stable, resistant to environmental changes and show no significant isomerization. ATTO 647N shows red fluorescence (λ_{abs} (max): 645nm, λ_{em} (max): 669nm, ϵ_{max} : 120'000).

UniProt ID

O00220

Worry-free Guarantee

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

Last modified: May 29, 2024



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