# 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-kB.

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 GenelD)

**Host** Mouse

Immunogen Recombinant human TRAIL-R1.

lgG1

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.

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 (λabs (max): 645nm, λem (max): 669nm, Emax: 120'000).

**UniProt ID** 

O00220

**Worry-free Guarantee** 

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-

uk@enzolifesciences.com