# DR3 (human):Fc (human), (recombinant)

The DR3 receptor (WSL-1, Apo-3, LARD, TRAMP, TR3) is a member of the tumor necrosis factor receptor (TNF-R) family. DR3 is a type I transmembrane receptor. It is 339 aa long and contains 4 extracellular cysteine-rich motifs in its extracellular domain, while its cytoplasmic region contains a death domain. DR3 is abundantly expressed on thymocytes and T and B lymphocytes. Overexpression of DR3 mediates activation of nuclear factor NF-κB and induces apoptosis. Besides the induction of apoptosis, DR3 has been reported to regulate negative selection during thymocyte development.

# **Ordering Information**

**Order Online** »

ALX-522-040-C050

50µg

Manuals, SDS & CofA

**View Online** »

### **Handling & Storage**

**Use/Stability** Stable for at least 6 months after receipt when stored at -20°C. After reconstitution, it is

recommended to make appropriate aliquots and to store them at -20°C.

**Handling** Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

# Regulatory Status RUO - Research Use Only

#### **Product Details**

Alternative Name WSL-1, Apo-3, LARD, TRAMP, TR3, TNFRSF 12,

TNFRSF 25, Death receptor 3

Application Notes Can be used together with anti-human IgG1 PAb to detect

TL1A.

Concentration 1mg/ml after reconstitution.

**Formulation** Lyophilized. Contains PBS.

**MW** ~55kDa (under reducing conditions).

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 cysteine-rich region of

human DR3 (aa 25-199) is fused to the Fc portion of

human IgG1.

**Specificity** Binds human TL1A.

UniProt ID Q93038



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

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

Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com