

# CD40L, soluble (human) ELISA kit (high sensitivity)

The CD40L, soluble (human) ELISA kit (high sensitivity) is a complete assay for the quantitative determination of soluble human CD40L (CD154) concentration. This kit provides an ultra-sensitive measurement of CD40L, detecting as little as 0.005ng/ml. This kit is to be used with cell culture supernatants, serum and other body fluids from human origin.

Citations: 6

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Handling** Store controls at -20°C. Store rest of the kit at +4°C.

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** CD154, TNFSF 5

**Application** ELISA

**Application Notes** For the quantitative determination of soluble human CD40L (CD154) in cell culture supernatants, serum and other body fluids.

**Assay Time** 3 hours + 10 minutes

**Compatibility** This product is compatible with the [Absorbance 96 Plate Reader](#).

**Quantity** 96 wells (~80 tests).

**Sensitivity** 0.005ng/ml (range 0.08 to 5ng/ml)

**Species Reactivity** Human

**UniProt ID** P29965

**Wavelength** 450 nm

**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-fr@enzolifesciences.com](mailto:info-fr@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)