

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 »](#)

Ordering Information

[Order Online »](#)

ALX-850-311-KI01	96 wells
------------------	----------

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



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)