

Product datasheet

Human TRAIL ELISA Kit (CD253) ab46074

[1 References](#) [3 Images](#)

Overview

Product name Human TRAIL ELISA Kit (CD253)

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
A	40	1500pg/ml	82.9	5.5%
B	33	116.3pg/ml	8.7	7.4%

Inter-assay

Sample	n	Mean	SD	CV%
A	12	1619pg/ml	62.8	3.8%
B	12	162.8pg/ml	13.1	8%

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity < 64 pg/ml

Range 93.75 pg/ml - 3000 pg/ml

Assay time 3h 45m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview

Abcam's TRAIL (CD253) Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of TRAIL in cell culture supernatants, buffered solutions or Human serum, plasma, and other body fluids.

A monoclonal antibody specific for TRAIL has been coated onto the wells of the microtiter strips provided. Samples, including standards of known TRAIL concentrations, control specimens or unknowns are pipetted into these wells. During the first incubation, the standards or samples and a biotinylated monoclonal antibody specific for TRAIL are simultaneously incubated. After washing, the enzyme Streptavidin-HRP, that binds the biotinylated antibody is added, incubated and washed. A TMB substrate solution is added which acts on the bound enzyme to induce a

colored reaction product. The intensity of this colored product is directly proportional to the concentration of TRAIL present in the samples.

This kit will recognize both endogenous and recombinant Human TRAIL.

Tested applications

Suitable for: Sandwich ELISA

Platform

Microplate

Properties

Storage instructions

Store at +4°C. Please refer to protocols.

Components	Identifier	1 x 96 tests	2 x 96 tests
10X Standard Diluent Buffer	Black	1 x 25ml	1 x 25ml
200X Wash Buffer	White	1 x 10ml	2 x 10ml
Biotinylated Antibody Diluent	Red	1 x 7ml	1 x 13ml
Biotinylated anti-TRAIL	Red	1 x 400µl	2 x 400µl
Chromogen TMB Substrate Solution		1 x 11ml	1 x 24ml
Control	Silver	2 vials	4 vials
HRP Diluent	Red	1 x 23ml	1 x 23ml
Plastic Plate Covers		2 units	4 units
Stop Reagent	Black	1 x 11ml	2 x 11ml
Streptavidin-HRP		2 x 5µl	4 x 5µl
TRAIL Microplate (12 x 8 well strips)		1 unit	2 units
TRAIL Standard (Lyophilized)	Yellow	2 vials	4 vials

Relevance

TRAIL is a cytokine that belongs to the tumor necrosis factor (TNF) ligand family. It preferentially induces apoptosis in transformed and tumor cells, but does not appear to kill normal cells although it is expressed at a significant level in most normal tissues. TRAIL binds to several members of TNF receptor superfamily including TNFRSF10A/TRAILR1, TNFRSF10B/TRAILR2, TNFRSF10C/TRAILR3, TNFRSF10D/TRAILR4, and possibly also to TNFRSF11B/OPG. The activity of TRAIL may be modulated by binding to the decoy receptors TNFRSF10C/TRAILR3, TNFRSF10D/TRAILR4, and TNFRSF11B/OPG that cannot induce apoptosis. The binding of TRAIL to its receptors has been shown to trigger the activation of MAPK8/JNK, caspase 8, and caspase 3.

Cellular localization

Cell Membrane; single-pass type II membrane protein.

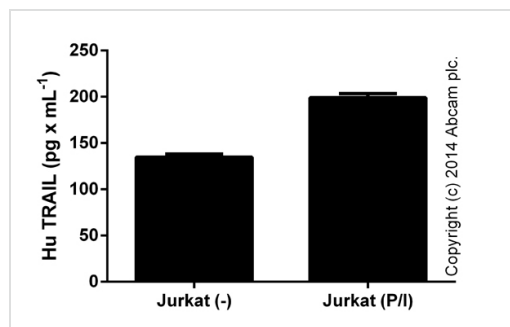
Applications

Our [Abpromise guarantee](#) covers the use of **ab46074** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

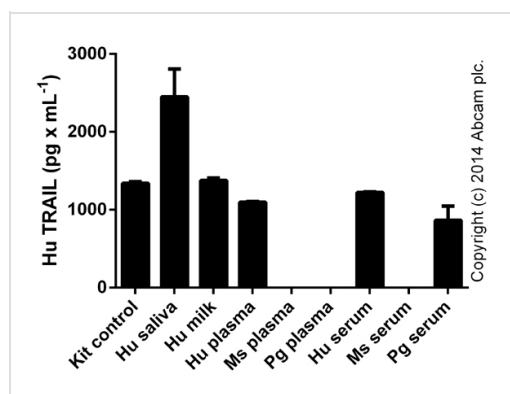
Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent concentration.

Images



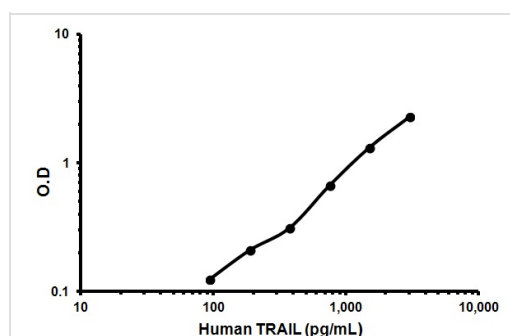
TRAIL measured in undiluted cell culture supernatants from control (-) or stimulated ((P/I); 48 hours with 50 ng x mL⁻¹ PMA ([ab120297](#)) and 1 uM ionomycin ([ab120116](#))) Jurkat cells (duplicates +/- SD).

ELISA: TRAIL (CD253) Human ELISA Kit (ab46074)



TRAIL measured in undiluted biological fluids (duplicates +/- SD).

ELISA: TRAIL (CD253) Human ELISA Kit (ab46074)



Representative Standard Curve using ab46074

Typical Standard Curve

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours

- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <http://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors