

## Product datasheet

# N-epsilon-(Carboxyethyl) Lysine (CEL) Assay Kit ab238540

[1 Image](#)

### Overview

<b>Product name</b>	N-epsilon-(Carboxyethyl) Lysine (CEL) Assay Kit
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Serum, Plasma, Cell Lysate, Purified protein
<b>Assay type</b>	Competitive
<b>Range</b>	0.1 µg/ml - 25 µg/ml
<b>Product overview</b>	N-epsilon-(Carboxyethyl) Lysine (CEL) Assay Kit (ab238540) is designed for the rapid detection and quantitation of CEL protein adducts.

First, a CEL conjugate is coated on the ELISA plate. The unknown CEL protein samples or CEL-BSA standards are then added to the CEL conjugate preabsorbed plate. After a brief incubation the anti-CEL monoclonal antibody is added, followed by an HRP conjugated secondary antibody. The content of CEL protein adducts in unknown samples is determined by comparison with the pre-determined CEL-BSA standard curve. Despite the structure similarity between CEL and CML, the anti-CEL specific antibody in this kit will not cross react with CML protein adducts.

**Platform** Microplate reader

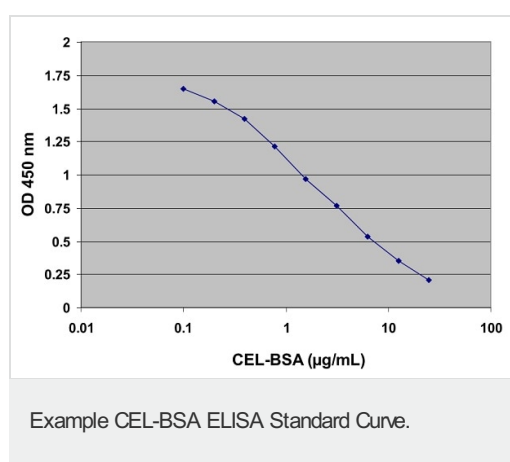
### Properties

**Storage instructions** Please refer to protocols.

Components	96 tests
100X Conjugate Diluent	1 x 300µl
10X Wash Buffer	1 x 100ml
Protein Binding Strip Well Plate	1 unit
Anti-CEL Antibody (1000X)	1 x 10µl
Assay Diluent	1 x 50ml

Components	96 tests
CEL Conjugate	1 x 20µl
CEL-BSA Standard	1 x 30µl
Secondary Antibody, HRP Conjugate (1000X)	1 x 20µl
Stop Solution	1 x 12ml
Substrate Solution	1 x 12ml

## Images



Typical CEL ELISA results.

This data is for reference only and should not be used to interpret actual results.

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

## 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 <https://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