

## Product datasheet

# Protein Carbonyl Content Assay Kit (Fluorometric) ab235631

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

### Overview

<b>Product name</b>	Protein Carbonyl Content Assay Kit (Fluorometric)
<b>Detection method</b>	Fluorescent
<b>Sample type</b>	Serum, Plasma, Cell Lysate
<b>Assay type</b>	Quantitative
<b>Sensitivity</b>	1 mg/ml
<b>Species reactivity</b>	<b>Reacts with:</b> Human
<b>Product overview</b>	Protein Carbonyl Content Assay Kit (Fluorometric) (ab235631) uses Fluorescein-5-Thiosemicarbazide (FTC), a fluorescent probe which covalently reacts with oxidized residues (i.e. Cysteine, Lysine, Arginine, Histidine and Aspartic Acid) on proteins.

Protein carbonyl content is determined by the generation of a stable fluorometric signal (Ex/Em 485/535 nm), which is compared to the protein concentration determined in a BCA Assay to quantitate nmoles of carbonyl/mg protein.

The fluorometric protein carbonyl content assay kit is simple, requires no harsh chemicals, can quantitate carbonyls in serum or plasma and produces more reliable and reproducible results than the comparable colorimetric assays. It can detect carbonyl groups in samples with protein concentrations as low as 1 mg/mL.

**Notes** This product is manufactured by BioVision, an Abcam company and was previously called K563 Protein Carbonyl Content Assay Kit (Fluorometric). K563-100 is the same size as the 100 test size of ab235631.

#### Related products

Other Protein Carbonyl Content assays include:

- [Protein Carbonyl Content Assay \(DNPH\) ab126287](#) (most popular protein carbonyl content assay)
- [Protein Carbonyl Content ELISA \(DNPH\) ab238536](#)
- [Protein Carbonyl Content Western Blot Assay ab178020](#)

Review the [oxidative stress marker and assay guide](#) to learn about more assays for oxidative stress.

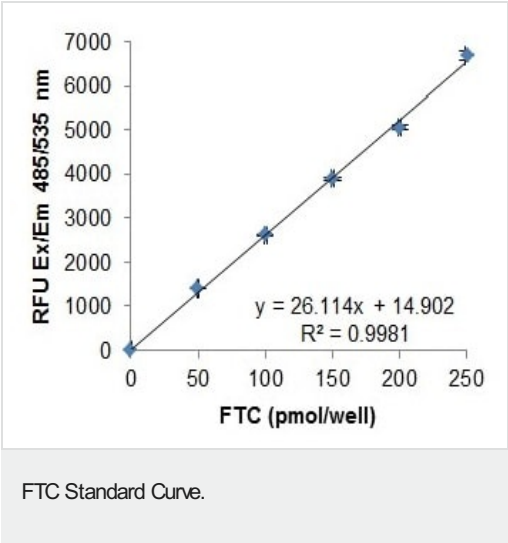
**Platform** Microplate reader

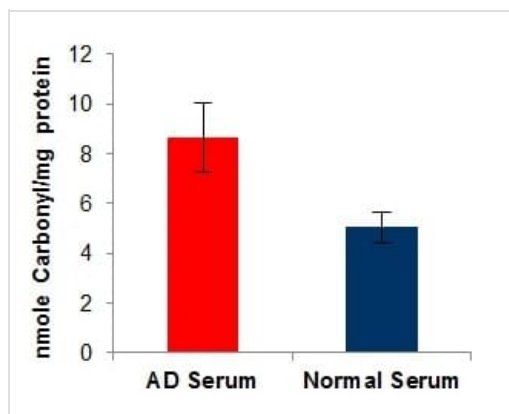
Properties

Storage instructions                      Store at -20°C. Please refer to protocols.

Components	100 tests
Protein Carbonyl Assay Buffer	1 x 20ml
100% TCA Solution	1 x 20ml
Streptozocin Solution	1 x 1ml
Guanidine Solution	1 x 20ml
Carbonyl Sample Dilution Buffer	1 x 50ml
Fluorescein-5-Thiosemicarbazide (FTC)	1 x 200µl
AGE-BSA Positive Control Protein	1 x 250µl

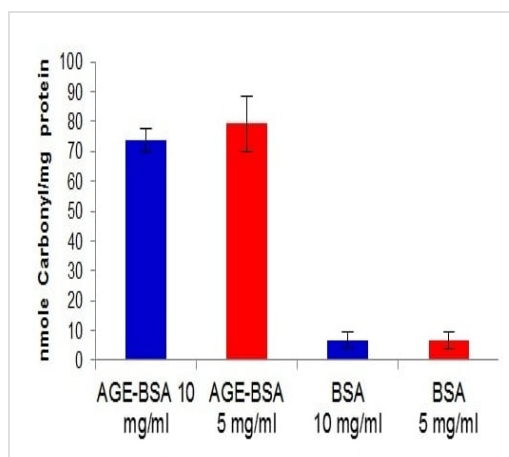
Images





50  $\mu$ L of human serum from patients with Alzheimer's Disease (AD) or healthy individuals were assayed following kit protocols.

Evaluation of carbonyl content in human serum.



Comparison of AGE-BSA with untreated BSA illustrating the 7-fold increase of carbonyls on AGEs as compared to BSA.

**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