

Product datasheet

Deproteinizing Sample Preparation Kit - TCA ab204708

★★★★★ [1 Abreviews](#) [57 References](#) [2 Images](#)

Overview

Product name	Deproteinizing Sample Preparation Kit - TCA
Sample type	Urine, Whole Blood, Tissue Extracts, Tissue Culture Media, Tissue Lysate
Assay time	0h 25m
Product overview	Deproteinizing Sample Preparation Kit - TCA (ab204708) offers a simple method to prepare samples for use in assays to measure small molecules, such as lipids, metabolites, metal ions etc.

This kit is useful where endogenous enzymes and other proteins would interfere with the assay, or would act directly to affect the level of the analyte to be measured, or would bind the analyte.

In this kit, trichloroacetic acid (TCA) is used to precipitate proteins by drastically lowering the sample pH. After removal of precipitated proteins, the pH of the sample is neutralized with neutralization buffer that is provided in the kit. Small molecules present in the sample can then be easily analyzed without interference from endogenous enzymes and other proteins .

The kit is easy to follow, convenient, and can be utilized for the preparation of a large number of samples in parallel. Samples prepared using this kit can be directly used in various bioassays.

Notes	This product is manufactured by BioVision, an Abcam company and was previously called K823 Deproteinizing Sample Preparation Kit II. K823-200 is the same size as the 200 test size of ab204708.
--------------	--

Tested applications	Suitable for: Purification
----------------------------	-----------------------------------

Properties

Storage instructions	Store at room temperature. Please refer to protocols.
-----------------------------	---

Components	200 tests
Neutralization Buffer I	1 x 4ml
Trichloroacetic Acid Solution	1 x 3ml

Applications

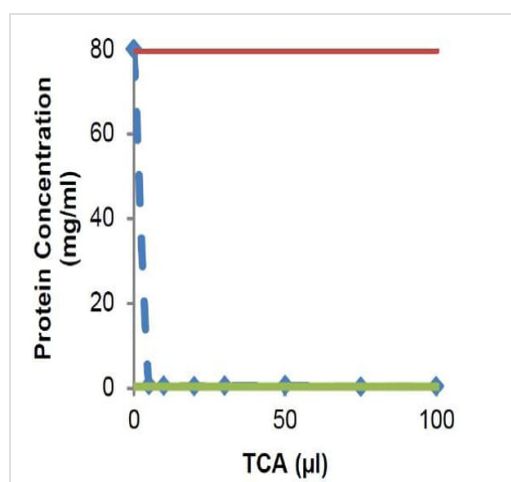
The Abpromise guarantee

Our **Abpromise guarantee** covers the use of ab204708 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
Purification		Use at an assay dependent concentration.

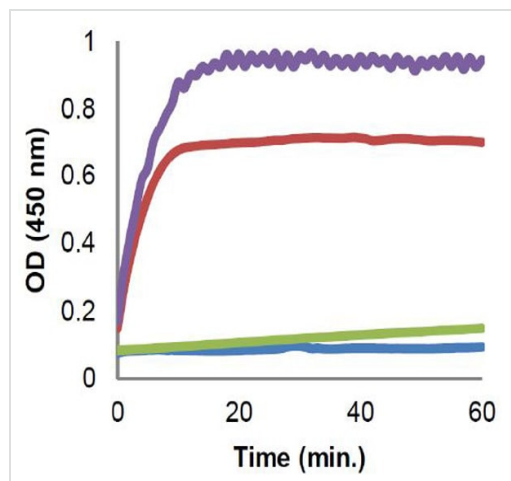
Images



ab 204708

Effect of TCA amount on deproteinization:

Aliquots of pooled human serum (off the clot, 100 µL) were treated with different amounts of TCA following the kit protocol. Estimated protein concentration before deproteinization: 7.9 g/dL. More than 99% of protein was removed using less than 10 µL of TCA/100 µL human serum. Red solid line = Total protein; blue dotted line = sample treated with TCA; green solid line = sample treated with spin column.



ab 204708

Comparison of analysis of lactate:

Human serum samples (100 µL) were analyzed for lactate concentration using L-Lactate Assay Kit (Colorimetric) ([ab65331](#)). Untreated samples (purple line) show higher signal when compared to TCA treated samples (red line). However, background signal is very similar in both untreated (green line) and treated (blue line) samples.

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