

Oxalate Assay Kit (Colorimetric) ab196990

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

Overview

Product name	Oxalate Assay Kit (Colorimetric)
Detection method	Colorimetric
Sample type	Urine, Serum, Plasma, Tissue
Assay type	Quantitative
Sensitivity	< 20 μ M
Assay time	2h 20m
Species reactivity	Reacts with: Mammals, Other species
Product overview	<p>Oxalate Assay kit (Colorimetric) (ab196990) is an easy-to-use, sensitive and high throughput adaptable kit. In this assay, Oxalate reacts with oxalate converter and oxalate enzyme mix to form an intermediate that reacts in turn with a highly specific probe to generate color that can be detected at OD=450 nm.</p> <p>This product can detect Oxalate levels lower than 20 μM.</p> <p>Oxalate assay protocol summary:</p> <ul style="list-style-type: none">- add samples and standards to wells- add oxalate convertor and incubate for 1 hr- add reaction mix and incubate for 1 hr- analyze with microplate reader
Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K663 Oxalate (Oxalic Acid) Colorimetric Assay Kit. K663-100 is the same size as the 100 test size of ab196990.</p> <p>Oxalate ($C_2O_4^{2-}$), in the form of Oxalic acid is present in many foods and beverages (e.g. spinach, tea etc.). It accumulates in many plant tissues and play role in regulating pH, osmosis and calcium storage. In animals, oxalate is either absorbed from dietary intake or produced from glycolate metabolism in liver. Under normal conditions, the daily oxalate load can be excreted by kidney. However, hereditary defects can cause an increased level of oxalate, which leads to hyperoxaluria and results in the formation of kidney stones. Therefore, measurement of oxalate level is useful for the prevention, diagnosis and monitoring of kidney stones.</p>
Platform	Microplate reader

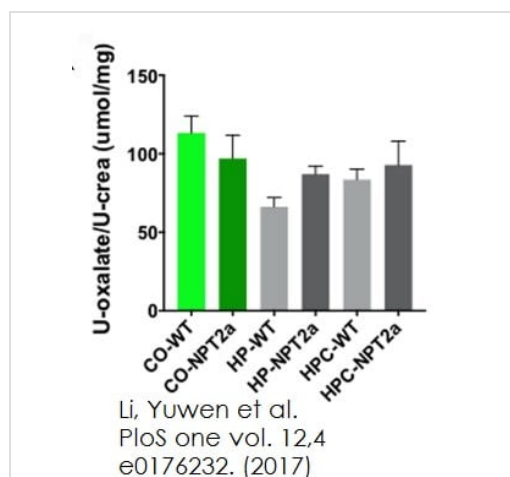
Properties

Storage instructions

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

Components	100 tests
Assay Buffer XXI	1 x 25ml
Developer Solution III	1 vial
Enzyme Mix IX	1 vial
Oxalate Converter	1 x 0.2ml
Buffer II	1 x 15ml
Oxalate Standard	1 vial

Images

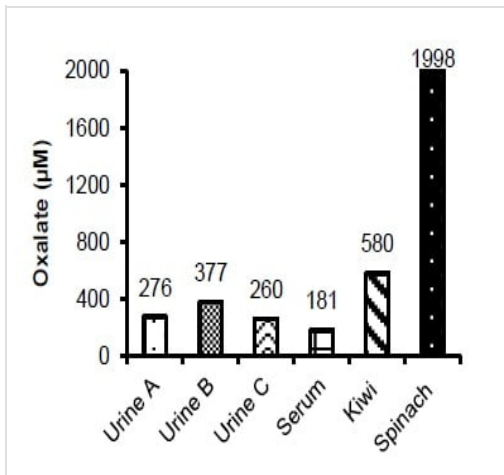


8 weeks old mice were placed for 10 weeks on special egg-white based diets: HPC diet (High phosphate and calcium diet; 20% lactate, 2% calcium, 1.25% phosphate); HP diet (High phosphate diet; 0.6% calcium, 1.20% phosphate); CO diet (Control diet; 0.6% calcium, 0.3% phosphate); WT: wild type; Npt2a: Npt2a^{-/-} mice. Oxalate was measured using 196990

Functional Studies - Oxalate Assay Kit

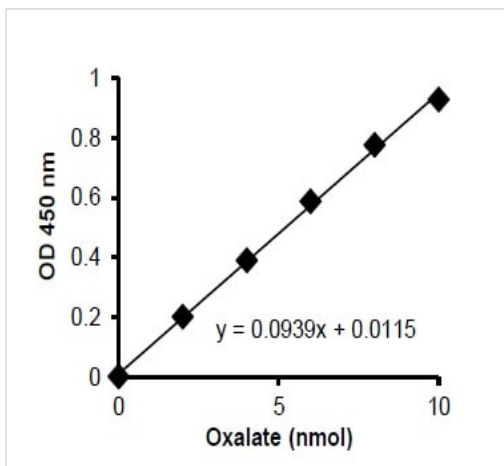
(Colorimetric) (ab196990)

Li, Yuwen et al., PloS one?vol. 12,4 e0176232., Fig 2,
doi:10.1371/journal.pone.0176232



Oxalate concentration in normal human urine (3 donors, 10 µL sample), normal human serum (10 µL), Kiwi lysate (10 µL) and Spinach lysate (10 µL).

Example data



Oxalate Standard Curve.

Example data

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