

# Glutamate Oxidase Assay Kit (Fluorometric - Near Infrared) ab138885

[1 References](#) [画像数 1](#)

### 製品の概要

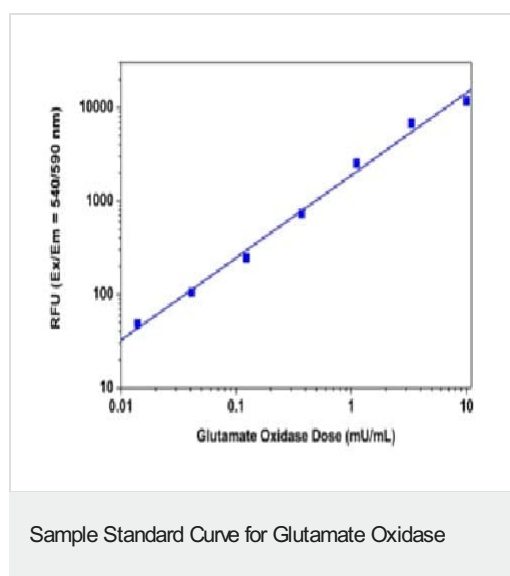
製品名	Glutamate Oxidase Assay Kit (Fluorometric - Near Infrared)
検出方法	Colorimetric/Fluorometric
サンプルの種類	Cell Lysate
アッセイタイプ	Enzyme activity
検出感度	= 40 uU/ml
種交差性	交差種: Mammals, Other species
製品の概要	<p>Glutamate Oxidase Assay Kit (Fluorometric - Near Infrared) (ab138885) provides a quick and an ultrasensitive method for the measurement of glutamate oxidase in solution and in cell lysates. In the assay, L-glutamic acid is oxidized to α-ketoglutarate, NH<sub>3</sub> and H<sub>2</sub>O<sub>2</sub> by glutamate oxidase.</p> <p>The kit uses our AbRed Indicator substrate which can react with H<sub>2</sub>O<sub>2</sub> when catalyzed by horseradish peroxidase (HRP) to generate the highly fluorescent product. The signal can be read with either a fluorescence microplate reader at Ex/Em = 530-570 nm/590 -600 nm (optimal Ex/Em = 540 nm/590 nm) or an absorbance microplate reader at 576±5 nm. With this product, we have detected as little as 40 μU/mL glutamate oxidase in a 100 μL reaction volume. It can be performed in a convenient 96-well or 384-well microtiter-plate format and easily adapted to automation without a separation step.</p>
特記事項	<p>Glutamate oxidase specifically catalyzes the oxidative deamination of L-glutamate in the presence of water and oxygen with the formation of ketoglutarate, ammonia, and hydrogen peroxide. Glutamate oxidase can be used as an analytic reagent and as a basis for developing biosensors for the determination of L-glutamate, L-glutamine, ammonia, and creatinine. These biosensors can be used in clinical biochemistry for the determination of glutamate-pyruvate transaminase and glutamate-oxalacetate transaminase in biological fluids, which makes the early diagnosis of heart and liver diseases possible.</p>
試験プラットフォーム	Microplate reader

### 製品の特性

保存方法 Store at -20°C. Please refer to protocols.

内容	200 tests
AbRed Indicator	1 vial
Assay Buffer	1 x 20ml
DMSO	1 x 200µl
Glutamate Oxidase Standard	1 vial
Glutamic Acid	1 x 3.4mg
Horseradish Peroxidase	1 vial

## 画像



Glutamate oxidase dose response was measured on a 96- well black plate with ab138885. As low as 40 µU/mL glutamate oxidase can be detected with 60 minutes incubation time (n=3).

**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.co.jp/abpromise> or contact our technical team.

## Terms and conditions

---

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