

# Arginase Activity Assay Kit (Colorimetric) ab180877

[19 References](#) [2 Images](#)

### Overview

<b>Product name</b>	Arginase Activity Assay Kit (Colorimetric)
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Tissue, Adherent cells, Suspension cells
<b>Assay type</b>	Enzyme activity (quantitative)
<b>Sensitivity</b>	< 0.2 U/ml
<b>Assay time</b>	1h 00m
<b>Species reactivity</b>	<b>Reacts with:</b> Mammals, Other species
<b>Product overview</b>	Arginase Activity Assay Kit (Colorimetric) ab180877 is a simple, sensitive and rapid assay to quantify arginase activity.

In the arginase assay protocol, arginase reacts with arginine and undergoes a series of reactions that form an intermediate that react stoichiometrically with the probe to generate a colored product that can be detected at OD = 570 nm.

The kit can detect less than 0.2 U/L Arginase activity in 96-well assay format.

Arginase assay protocol summary:

- add samples and standards to wells
- add arginase substrate mix to sample wells only
- incubate for 20 min
- add arginase assay reaction mix to all wells
- analyze with microplate reader in kinetic mode for 10-30 min

**Notes** This product is manufactured by BioVision, an Abcam company and was previously called K755 Arginase Activity Colorimetric Assay Kit. K755-100 is the same size as the 100 test size of ab180877.

Arginase (EC 3.5.3.1) is the final enzyme of the Urea Cycle. It converts L-arginine into urea and L-ornithine and plays an important role in removing ammonium ion from the body.

**Platform** Microplate reader

### Properties

**Storage instructions**

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

Components	100 tests
Arginase Positive Control	1 vial
Arginine	1 vial
Assay Buffer XIV	1 x 25ml
Converter Enzyme V	1 vial
Developer I	1 vial
Development Enzyme Mix I	1 vial
H2O2 Standard	1 x 100µl
OxiRed Probe	1 x 200ml

**Pathway**

Nitrogen metabolism; urea cycle; L-ornithine and urea from L-arginine: step 1/1.

**Involvement in disease**

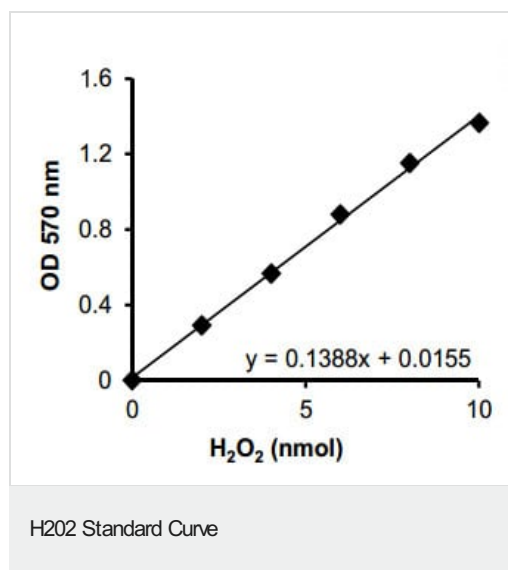
Defects in ARG1 are the cause of argininemia (ARGIN) [MIM:207800]; also known as hyperargininemia. Argininemia is a rare autosomal recessive disorder of the urea cycle. Arginine is elevated in the blood and cerebrospinal fluid, and periodic hyperammonemia occurs. Clinical manifestations include developmental delay, seizures, mental retardation, hypotonia, ataxia, progressive spastic quadriplegia.

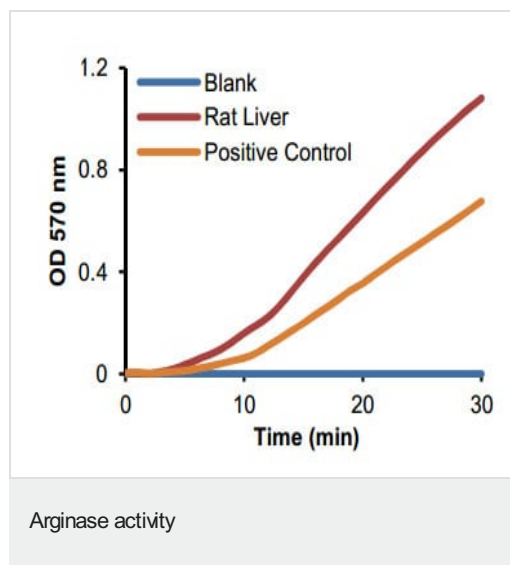
**Sequence similarities**

Belongs to the arginase family.

**Cellular localization**

Cytoplasm.

**Images**H<sub>2</sub>O<sub>2</sub> Standard Curve. Assay performed following kit protocol.



Arginase activity in rat liver lysate (3 µg) & Positive Control (2 µL).  
Assays were performed following the kit protocol.

**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