

## Product datasheet

# Gluconokinase Activity Assay Kit (Colorimetric) ab211092

[3 Images](#)

### Overview

<b>Product name</b>	Gluconokinase Activity Assay Kit (Colorimetric)
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Cell Lysate, Tissue Lysate
<b>Assay type</b>	Enzyme activity (quantitative)
<b>Sensitivity</b>	< 100 µU
<b>Species reactivity</b>	<b>Reacts with:</b> Mammals, Other species
<b>Product overview</b>	Gluconokinase Activity Assay Kit (Colorimetric) (ab211092) provides a simple, sensitive and convenient method for detecting gluconokinase activity in a variety of samples such as cell and tissue lysates, and even in prokaryotes such as <i>E. coli</i> . In this assay, Gluconokinase converts Gluconate into 6-Phosphate-D-Gluconate in an ATP dependent manner. 6-Phosphate-D-Gluconate and ADP in turn undergo a series of reactions to form an intermediate, which reacts with the probe to form a colored product with strong absorbance (OD 450 nm). The detection limit of this assay < 0.1 mU.

<b>Notes</b>	<p>This product is manufactured by BioVision, an Abcam company and was previously called K319 Gluconokinase (GntK) Activity Assay Kit (Colorimetric). K319-100 is the same size as the 100 test size of ab211092.</p> <p>Gluconokinase (GntK, ATP:D-gluconate 6-phosphotransferase, Gluconate Kinase; EC:2.7.1.12) is a key enzyme for degradation of gluconate. In <i>E. coli</i> and yeast, Gluconokinase phosphorylates gluconate into 6-Phosphate-D-Gluconate in an ATP dependent manner. Via the Hexose Monophosphate Shunt (HMS) pathway, 6- Phosphate-D-Gluconate generates ribose-5-phosphate, which is critical for nucleotides and nucleic acid synthesis.</p> <p>Despite the widespread use of D-gluconate as acidity regulator in food and medicine, little is known of the mechanism of gluconate metabolism in humans.</p>
<b>Platform</b>	Microplate reader

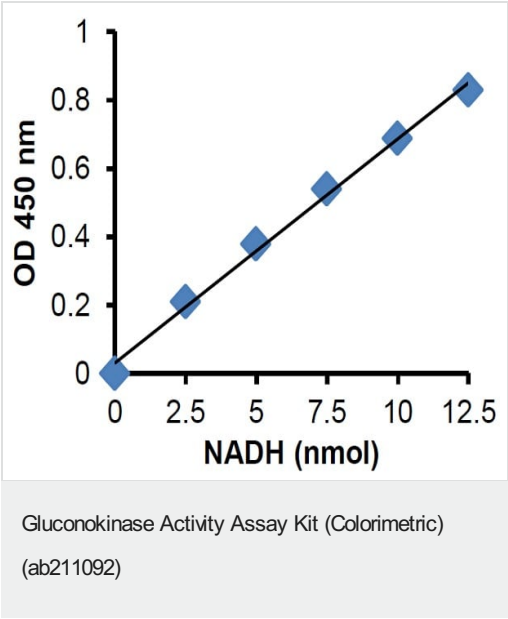
### Properties

<b>Storage instructions</b>	Store at -20°C. Please refer to protocols.
-----------------------------	--

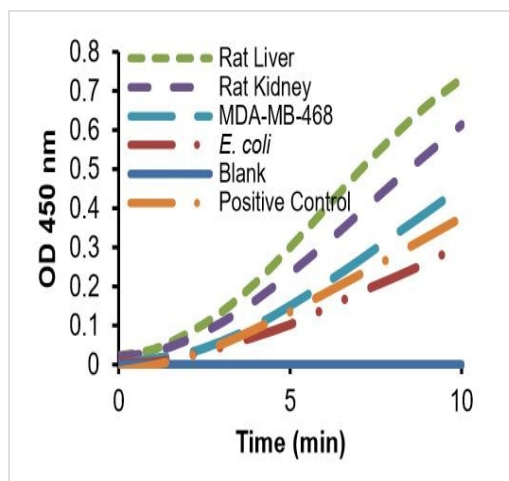
Components	100 tests
GntK Assay Buffer	1 x 25ml
ATP (40 $\mu$ mole)	1 vial
GntK Converting Enzyme (10 U)	1 vial
GntK Developer (10 U)	1 vial
GntK Positive Control (200 mU)	1 vial
GntK Probe (30 mg)	1 vial
GntK Substrate (50 $\mu$ mole)	1 vial
NADH Standard (0.5 $\mu$ mole)	1 vial

Pathway	Carbohydrate acid metabolism; D-gluconate degradation.
Sequence similarities	Belongs to the gluconokinase gntK/gntV family.

Images

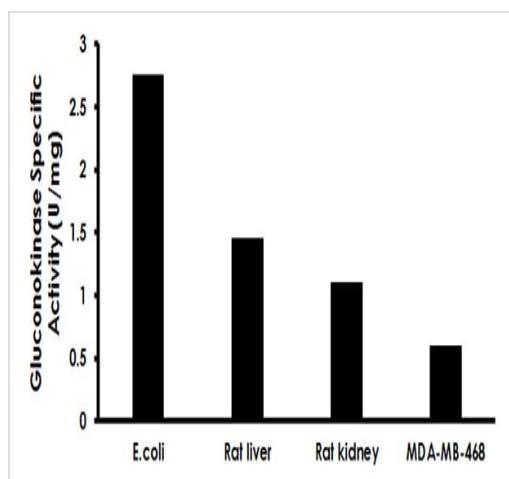


Typical NADH standard calibration curve.



Kinetic curves showing Gluconokinase activity detection in positive control (1 µL; included in kit), lysates from rat liver and kidney (1 µg, respectively), lysates from MDA-MB-468 cells (1 µg) and *E. coli* (0.1 µg).

Gluconokinase Activity Assay Kit (Colorimetric)  
(ab211092)



Gluconokinase specific activity in lysates from rat liver and kidney (1 µg, respectively), lysates from MDA-MB-468 cells (1 µg) and *E. coli* (0.1 µg).

Gluconokinase Activity Assay Kit (Colorimetric)  
(ab211092)

**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