



# Invertase Assay Kit

Catalog Number KA1629

100 assays

Version: 04

Intended for research use only

[www.abnova.com](http://www.abnova.com)

## Table of Contents

<b>Introduction .....</b>	<b>3</b>
Intended Use .....	3
Background .....	3
Principle of the Assay .....	3
<b>General Information .....</b>	<b>4</b>
Materials Supplied .....	4
Storage Instruction .....	4
Materials Required but Not Supplied .....	4
Precautions for Use .....	4
<b>Assay Protocol .....</b>	<b>5</b>
Assay Procedure .....	5
<b>Data Analysis.....</b>	<b>6</b>
Calculation of Results.....	6
<b>Resources.....</b>	<b>7</b>
References .....	7

## Introduction

### Intended Use

- ✓ Application
  - Invertase and sucrase activity determination in biological and environmental (e.g. soil) samples.
  - Evaluation and screening for invertase inhibitors.
- ✓ Features:
  - Safe: Non-radioactive assay.
  - Sensitive and accurate: As low as 0.007 U/L invertase activity can be quantified.
  - Homogeneous and convenient: "Mix-incubate-measure" type assay. No wash and reagent transfer steps are involved.
  - Robust and amenable to HTS: can be readily automated on HTS liquid handling systems for processing thousands of samples per day.

### Background

INVERTASE ( $\beta$ -fructofuranosidase, EC 3.2.1.26) is an enzyme that catalyzes the hydrolysis of sucrose to fructose and glucose. Invertases cleave at the O-C(fructose) bond, whereas a related enzyme sucrase (EC 3.2.1.48) cleaves at the O-C(glucose) bond. A wide range of microorganisms produce invertase and can, thus, utilize sucrose as a nutrient. Invertase assay finds wide applications in environmental (e.g. soil), agricultural and food (confectionery) industry.

### Principle of the Assay

The Invertase Assay Kit provides a convenient and ultra-sensitive colorimetric and fluorimetric means to measure invertase activity. In the assay, invertase cleaves sucrose, resulting in the formation of fructose and glucose, which is determined by a colorimetric (570 nm) or fluorimetric method ( $\lambda_{em/ex}$  = 585/530 nm). The assay is simple, sensitive, stable and high-throughput adaptable.

## General Information

### Materials Supplied

List of component

Component	Amount
10x Reaction Buffer: (pH 4.5)	12 mL
Assay Buffer	10 mL
Glucose Standard	1 mL
10x Sucrose	1.5 mL
Enzyme Mix	120 µL
Dye Reagent	120 µL

### Storage Instruction

Store all components at -20°C. Shelf life of six months after receipt.

### Materials Required but Not Supplied

- ✓ Pipetting devices
- ✓ Centrifuge tubes
- ✓ Clear or black flat bottom 96-well plate (e.g. Corning Costar)

### Precautions for Use

- ✓ Precautions

Reagents are for research use only. Normal precautions for laboratory reagents should be exercised while using the reagents.

## Assay Protocol

### Assay Procedure

Interference: thiols ( $\beta$ -mercaptoethanol, dithioerythritol etc) at  $> 10 \mu\text{M}$  interfere with this assay and should be avoided. Glucose, if present in the sample, should be removed by dialysis or membrane filtration.

#### 1. Assay Preparation

Prior to assay, bring all components to room temperature, briefly centrifuge tubes before opening. Dilute the provided 10x Reaction Buffer and 10x Sucrose to 1-fold by mixing 1 vol of the reagent with 9 vol of  $\text{dH}_2\text{O}$ . Use the diluted reagents for all assays.

For glucose standard curve, mix 5  $\mu\text{L}$  Glucose Standard with 828  $\mu\text{L}$   $\text{dH}_2\text{O}$  (final 100  $\mu\text{M}$ ). Dilute as follows and transfer 40  $\mu\text{L}$  standards to separate wells in a clear flat-bottom 96-well plate.

No	100 $\mu\text{M}$ Std + $\text{H}_2\text{O}$	Vol ( $\mu\text{L}$ )	Glucose ( $\mu\text{M}$ )
1	100 $\mu\text{L}$ + 0 $\mu\text{L}$	100	100
2	60 $\mu\text{L}$ + 40 $\mu\text{L}$	100	60
3	30 $\mu\text{L}$ + 70 $\mu\text{L}$	100	30
4	0 $\mu\text{L}$ + 100 $\mu\text{L}$	100	0

Sample: transfer 40  $\mu\text{L}$  sample to separate wells of the plate. As a sample control, use 40  $\mu\text{L}$  diluted Reaction Buffer.

#### 2. Enzyme Reaction

Add 5  $\mu\text{L}$  of the diluted Sucrose to each well. Tap plate to mix. Incubate 20 min at desired temperature (e.g. 30°C).

#### 3. Glucose Determination

Prepare enough Working Reagent in bulk.

For each well, mix 95  $\mu\text{L}$  Assay Buffer, 1  $\mu\text{L}$  Enzyme Mix, 1  $\mu\text{L}$  Dye Reagent. Add 90  $\mu\text{L}$  Working Reagent to each well. Immediately tap plate to mix.

Incubate for 20 min in the dark. Read OD570nm.

*Note: the procedure for fluorimetric assays is the same except that (1) a black flat-bottom 96-well plate is used, (2) glucose standards should be at 20, 12, 6 and 0  $\mu\text{M}$  and that fluorescence intensity at  $\lambda_{\text{em/ex}} = 585/530\text{nm}$  is measured.*

## Data Analysis

### Calculation of Results

Plot glucose standard curve and determine its Slope ( $\mu\text{M}^{-1}$ ). Invertase enzyme activity in the sample is calculated as

$$\text{Invertase Activity} = \frac{R_{\text{SAMPLE}} - R_{\text{CONTROL}}}{\text{Slope} \times t} \text{ (U/L)}$$

where  $R_{\text{SAMPLE}}$  and  $R_{\text{CONTROL}}$  are the OD or fluorescence values of the sample and sample control (i.e. Reaction Buffer).  $t$  is the incubation time (20 min).

Unit definition: one unit of invertase catalyzes the formation of 1  $\mu\text{mole}$  glucose per min at pH 4.5 under the assay conditions.

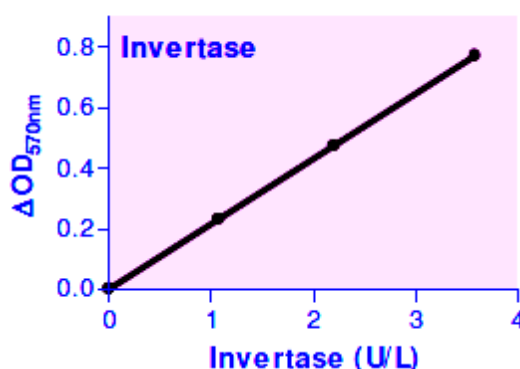
*Note: if the OD or fluorescence intensity is higher than the value for 100  $\mu\text{M}$  glucose (colorimetric assay) or 20  $\mu\text{M}$  (fluorimetric assay), dilute sample in 1-fold Reaction Buffer and repeat the assay. Multiply the result by the enzyme dilution factor.*

#### ✓ INVERTASE ASSAY IN SOIL SAMPLES

Soil samples can be directly assayed as follows. Weigh about 100 mg soil into a 1.5 mL Eppendorf tube. Add 880  $\mu\text{L}$  diluted Reaction Buffer and 120  $\mu\text{L}$  diluted sucrose. Mix thoroughly by homogenization and/or vortexing. Immediately remove 200  $\mu\text{L}$  mixture into a clean tube and centrifuge for 2 min at 14,000 rpm. Transfer 100  $\mu\text{L}$  clear supernatant into another clean tube and immediately freeze at  $-20^{\circ}\text{C}$ . This "time zero" sample serves as a sample control.

Incubate the invertase reaction for 1 hour at 30 or  $37^{\circ}\text{C}$  (Step 2). Centrifuge for 2 min at 14,000 rpm. Transfer 40  $\mu\text{L}$  clear supernatant and the above sample control for glucose determination (Step 3).

Example 1: purified yeast invertase



Example 2:

A 100 mg soil sample was assayed according to the above procedure. At the end of 1 hour enzyme reaction at  $30^{\circ}\text{C}$ , 58.4  $\mu\text{M}$  glucose was determined, which corresponds to an invertase activity of  $58.4 \mu\text{moles/L} \div 60 \text{ min} = 0.97 \text{ U/L}$ , or  $58.4 \mu\text{moles/L} \div (100 \text{ g/L} \times 1 \text{ hour}) = 0.58 \mu\text{moles} \times \text{g}^{-1} \times \text{hr}^{-1}$  or  $105.2 \mu\text{g glucose} \times \text{g}^{-1} \times \text{hr}^{-1}$ .

## Resources

### References

1. Huang, Y.H., Picha, D.H. and Johnson, C.E.(1998). An alternative method for enzymatic assay of plant invertases. J. Agric. Food Chem. 46 (8): 3158–3161.
2. Ewing, E.E. and McAdoo, M.H. (1971). An examination of methods used to assay potato tuber invertase and its naturally occurring inhibitor. Plant Physiol. 48: 366-370.
3. Silveira, M.C., Carvajal, E. and Bon, E.P. (1996). Assay for in vivo yeast invertase activity using NaF. Anal Biochem. 238: 26-28.