

## EnzyChrom™ Lactose Assay Kit (ELAC-100)

### Quantitative Colorimetric Lactose Determination

#### DESCRIPTION

Lactose (C<sub>12</sub>H<sub>22</sub>O<sub>11</sub>), also called milk sugar, is a disaccharide that consists of β-D-galactose and α/β-D-glucose through a β1-4 glycosidic linkage. Lactose is the major sugar and makes up 2–8% of milk. Simple, direct and high-throughput assays for lactose determination find wide applications. BioAssay Systems' assay uses specific enzyme-coupled reactions in which lactose is cleaved and the resulting galactose forms a colored product. The color intensity at 570nm or fluorescence intensity at 530nm/585nm is directly proportional to the lactose concentration in the sample.

#### KEY FEATURES

Use as little as 20 μL samples. Linear detection range in 96-well plate: 17 to 2000 μM lactose for colorimetric assays and 6 to 100 μM for fluorimetric assays.

#### APPLICATIONS

**Assays** of lactose in milk and other biological samples.

**Drug Discovery/Pharmacology:** effects of drugs on lactose metabolism.

**Food and Beverages:** lactose in food and beverages products.

#### KIT CONTENTS

<b>Assay Buffer:</b>	10 mL	<b>Enzyme Mix:</b>	Dried
<b>Dye Reagent:</b>	120 μL	<b>Lactase:</b>	Dried
<b>Standard (20 mM Lactose):</b> 1 mL			

**Storage conditions.** The kit is shipped on dry ice. Store all components at -20°C. Shelf life of 6 months after receipt.

**Precautions:** reagents are for research use only. Normal precautions for laboratory reagents should be exercised while using the reagents. Please refer to Material Safety Data Sheet for detailed information.

#### COLORIMETRIC PROCEDURE

*Note: (1) glycerol and SH-containing reagents (e.g. β-mercaptoethanol, dithiothreitol) are known to interfere in this assay and should be avoided in sample preparation. (2) For samples containing galactose, a sample blank is necessary (see Procedure); (3) This assay is based on a kinetic reaction. To ensure identical incubation time, addition of Working Reagent to standard and samples should be quick and mixing should be brief but thorough. Use of a multi-channel pipettor is recommended.*

**Sample treatment:** Milk samples should be cleared by mixing 600 μL milk with 100 μL 6 N HCl. Centrifuge 5 min at 14,000 rpm. Transfer 300 μL supernatant into a clean tube and neutralize with 50 μL 6 N NaOH. The neutralized supernatant is ready for assay (dilution factor  $n = 1.36$ ).

1. Equilibrate all components to room temperature. Reconstitute the Lactase and Enzyme mix with 120 μL dH<sub>2</sub>O. Reconstituted Lactase and Enzyme mix are stable for 3 months if stored at -20°C. During experiment, keep reconstituted Lactase and Enzyme Mix in a refrigerator or on ice.

2. **Standards and samples:** prepare 400 μL 2000 μM Standard by mixing 40 μL 20 mM standard with 360 μL dH<sub>2</sub>O. Dilute standard in dH<sub>2</sub>O as follows.

No	2000 μM STD + H <sub>2</sub> O	Vol (μL)	Lactose (μM)
1	100 μL + 0 μL	100	2000
2	80 μL + 20 μL	100	1600
3	60 μL + 40 μL	100	1200
4	40 μL + 60 μL	100	800
5	30 μL + 70 μL	100	600
6	20 μL + 80 μL	100	400
7	10 μL + 90 μL	100	200
8	0 μL + 100 μL	100	0

Transfer 20 μL standards and 20 μL samples into separate wells of a clear flat-bottom 96-well plate. *Note: if a sample is known to contain galactose, transfer 20 μL sample in duplicate (one sample and one sample blank).*

3. **Reaction.** For each reaction well, mix 85 μL Assay Buffer, 1 μL Lactase, 1 μL Enzyme Mix (vortex briefly before pipetting), and 1 μL Dye Reagent in

a clean tube. (Note: for the sample blanks, prepare a control Working Reagent which is the same except WITHOUT the 1 μL Lactase). Transfer 80 μL Working Reagent into each reaction (and control) well. Tap plate to mix. Incubate 30 min at room temperature.

4. Read optical density at 570nm (550-585nm).

#### FLUORIMETRIC PROCEDURE

For fluorimetric assays, the linear detection range is 6 to 100 μM lactose. Prepare 100 μM lactose standard by mixing 5 μL 20 mM standard with 995 μL H<sub>2</sub>O. Then dilute standards in H<sub>2</sub>O (see *Colorimetric Procedure*) to 100, 80, 60, 40, 30, 20, 10 and 0 μM.

1. Transfer 20 μL standards and 20 μL samples into separate wells of a black 96-well plate. Prepare Sample Blank if necessary.

2. Add 80 μL Working Reagent, tap plate to mix. Incubate 30 min.

3. Read fluorescence at  $\lambda_{ex} = 530\text{nm}$  and  $\lambda_{em} = 585\text{nm}$ .

**Notes:** If the calculated lactose concentration of a sample is higher than 2000 μM in colorimetric assay or 100 μM in fluorimetric assay, dilute sample in water and repeat the assay. Multiply result by the dilution factor  $n$ .

#### CALCULATION

Subtract blank value (water, #8) from the standard values and plot the  $\Delta OD$  or  $\Delta RFU$  against standard concentrations. Determine the slope and calculate the lactose concentration of Sample,

$$\text{Colorimetry: } [\text{Lactose}] = \frac{OD_{\text{SAMPLE}} - OD_{\text{BLANK}}}{\text{Slope}} \times n \quad (\mu\text{M})$$

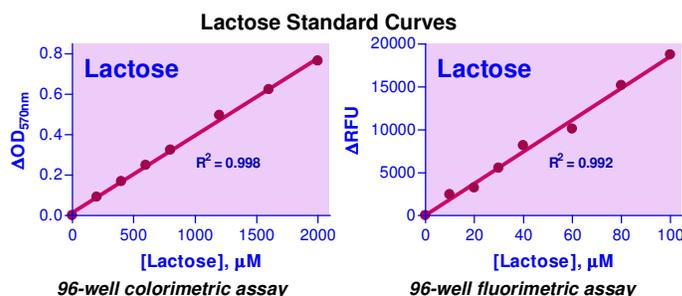
$$\text{Fluorimetry: } [\text{Lactose}] = \frac{RFU_{\text{SAMPLE}} - RFU_{\text{BLANK}}}{\text{Slope}} \times n \quad (\mu\text{M})$$

$OD_{\text{SAMPLE}}$ ,  $OD_{\text{BLANK}}$ ,  $RFU_{\text{SAMPLE}}$ ,  $RFU_{\text{BLANK}}$  are optical density and fluorescence values of the Sample and Blank. The Blank is water if there is no galactose, and Sample Blank if sample contains galactose.  $n$  is the dilution factor.

**Conversions:** 1 mM lactose equals 34.2 mg/dL, 0.0342% or 342 ppm.

#### MATERIALS REQUIRED, BUT NOT PROVIDED

Pipetting devices, centrifuge tubes, clear flat-bottom 96-well plates, optical density plate reader; black 96-well plates and fluorescence plate reader.



#### PUBLICATIONS

1. Xue, H et al. (2020). Lactose-induced chronic diarrhea results from abnormal luminal microbial fermentation and disorder of ion transport in the colon. *Frontiers in Physiology*, 11, 877.

2. Snyder, N. A., Palmer, M. V., Reinhardt, T. A., & Cunningham, K. W. (2019). Milk biosynthesis requires the Golgi cation exchanger TMEM165. *Journal of Biological Chemistry*, 294(9), 3181-3191.

3. Vabbilisetty, P and S Xue-Long (2014). Liposome surface functionalization based on different anchoring lipids via Staudinger ligation. *Organic and Biomolecular Chemistry* 12(8): 1237-44.

