

Cellometer[®]

Product Number: CS0-0110-500mL

Description: Cellometer ViaStain™ Yeast Dilution Buffer

Instrument (s): Cellometer X2, Vision10x, Spectrum 10x

Instruction Booklet: ViaStain™ Yeast Dilution Buffer



This product is for RESEARCH USE ONLY and is not approved for diagnostic or therapeutic use.

8001308 Rev. C

Left blank intentionally

Product

Part Number: **CS0-0110-500mL**

Description: Cellometer ViaStain™ Yeast Dilution Buffer

Description

The ViaStain™ Yeast Dilution Buffer in combination with the ViaStain™ Yeast Live / Dead Cell staining solution enables the user to quantitatively distinguish live and dead yeast in pure cultures and in cultures containing debris such as corn mash using the Cellometer system. The Dilution Buffer is used to dilute the yeast sample in order to promote and optimize live / dead fluorescent staining of the yeast cells.

Materials

Materials Supplied

1. Cellometer ViaStain™ Yeast Dilution Buffer (CS0-0110-500mL)*

*One 500 mL bottle of Yeast Dilution Buffer

Materials Required

1. Micro centrifuge tube
2. Pipettor
3. Cellometer counting chamber (SD100 or PD100)
4. Cellometer Vision10x (with Fluorescence Optical Module F101, VB-535-401, or equivalent and F304, VB-660-501, or equivalent), Spectrum 10x (with Fluorescence Optical Module S1-534-470), or a Cellometer X2

Procedure

1. Dilute corn mash sample 10-fold (by weight) in water and mix well.
2. Allow corn mash debris in diluted sample to settle for 30 sec.
3. Pipette 10 µL of diluted corn mash (while avoiding large debris in the mixture) and dispense into a microcentrifuge tube.
4. Add 10 µL of yeast dilution buffer to microcentrifuge tube and mix well by pipetting up and down 3 times.
5. Add 20 µL ViaStain™ Yeast Live / Dead Cell staining solution to microcentrifuge tube and mix well by pipetting up and down 3 times.
6. Incubate the stained corn sample for 1-2 min at room temperature.
7. Pipette sample up and down 3 times to mix and then load 20 µL into a counting chamber (if using SD100 slides, peel plastic film off both sides before loading).
8. Place loaded slide on a Kimwipe® and wait 1 min before inserting sample into instrument to allow sample to settle in the chamber.
9. Select the appropriate assay type for yeast viability measurement.
10. Preview bright field and fluorescent images.
11. Count.

Storage and Handling

Store the Yeast Dilution Buffer at 4°C. The ViaStain™ Yeast Dilution Buffer is a buffer at high pH; safety precautions must be taken when handling the solution. Please consult the Safety Data Sheet for more safety information.

Warranty

This product is for RESEARCH USE ONLY and is not approved for diagnostic or therapeutic use. Product is warranted to meet the specifications outlined in the Certificate of Analysis when stored and used according to the manufacturer's instructions. No other warranty, expressed or implied (such as merchantability, fitness for a particular purpose, or non-infringement) is granted. Warranty is valid until the expiration date stated on the product label. If no expiration is listed, the warranty is valid for 12 months from the date of product receipt.

Warranty will be void if product is stored incorrectly, the recommended protocol is not followed, or the product is used for a different application.

Ordering Information

When ordering with a Purchase Order:

Fax a copy of the order to 978-327-5341

Email a copy of the order to sales@nexcelom.com

When ordering with a Credit Card:

Visit www.shop.nexcelom.com and place your order