

Cellometer[®]

Product Number:

Full Product: CS1-0109-5mL

Sample Product: CS1-0109-S (Not for re-sale)

Description: ViaStain[™] PI Staining Solution

Instrument (s): Auto2000, X4, X1, X2, K2, Vision, Vision CBA, Spectrum

Product Insert: ViaStain[™] PI Staining Solution



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

8001306 Rev. D

Page Left Blank Intentionally

Description

The ViaStain™ PI Staining Solution enables the user to quantitatively distinguish live and dead cells in pure cultures using the Cellometer system. Propidium iodide (PI) is a membrane exclusion nucleic acid dye that only enters cells with compromised membranes (dead cells). When cell counting is performed using both bright-field and fluorescent imaging, cells that fluoresce in red are counted as dead, while the bright-field image provides the total cell count. The Cellometer software utilizes the total cell count and dead cell count to calculate percent viability.

Materials

Materials Supplied

1. CS1-0109-5mL (250 tests)
2. CS1-0109-S (50 Tests)

Materials Required

1. Micro centrifuge tube
2. Pipette
3. Cellometer counting chamber (SD100 or PD100)
4. Cellometer Auto 2000, Cellometer Auto X4, with Fluorescence Optical Module XB-595-501, Cellometer K2, Cellometer X1, Cellometer X2, Spectrum (with Fluorescent Optics Module S1-605-527 or S1-655-527) , or Cellometer Vision / Vision CBA, with Fluorescence Optical Module VB-595-502, VB-660-502 or equivalent

Procedure

1. Pipette 20 µL of cell sample into a micro centrifuge tube
2. Add 20 µL of PI staining solution to micro centrifuge tube and mix well by pipetting up and down at least 3 times.
3. Load 20 µL into a counting chamber (if using SD100 slides, peel plastic film off both sides before loading).
4. Insert loaded slide into the instrument.
5. Select the appropriate assay type for PI cell count and viability measurement.
6. Preview bright-field and fluorescent images.
7. Focus if necessary.
8. Count.

Storage and Handling

Store the PI Staining Solution at 2-8°C protected from light. AVOID FREEZING. Safety precautions must be taken when handling the solution. Please consult the Safety Data Sheet for more safety information, found on www.nexcelom.com/Products.

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 (for orders shipping in the United States):

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 (For orders shipping to a destination outside of the United States):

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