

Description

Part Number: **CCBM-011-2mL**
Description: Cellometer Check[™] Validation Bead Solution
Catalog Number: **CCBM-011-2mL**
Manufacturer Lot Number: **240401-01**

Characteristics

Color: Cellometer Green and Cellometer Red
Antimicrobial: **2mM NaN3**
Buffer: DI water + 0.003% Tween 20

Excitation/Emission Ranges

Full width half maximum
Cellometer Green Cellometer Red
Excitation: 475 - 515 Excitation: 460 - 560
Emission: 505 - 540 Emission: 540 - 660

Quality Control

Manufactured and tested according to SOP #: 8001826
Concentration:
Average concentration: 2.37×10^6 Beads/mL
Bead solution range = 2.01×10^6 to 2.73×10^6 Beads/mL
Viability:
% Viability = $\frac{\text{Live Beads}}{\text{Live Beads} + \text{Dead Beads}} = 76.13\%$
Specification: % Green FL Beads = 67.5 % to 82.5 %



Matthew Bularzik, Quality Engineer

04/01/2024

Recommended Storage Conditions:

Store the Validation Bead Solution at 2-8°C protected from light. **AVOID FREEZING!**

For research use only. Not intended for *in vivo* use.

If you have any further questions about this Certificate of Analysis, please contact us.
Nexcelom Bioscience | 360 Merrimack Street, Bldg 9 | Lawrence, MA 01843
T: 978.327.5340 | F: 978.327.5341 | E: info@nexcelom.com | www.nexcelom.com
8001344 Rev. C