

Cellometer[®] Check

Calibration Beads

Cat #	Bead diameter (μm)	Volume (mL)
CCBM-011-2mL	10	2

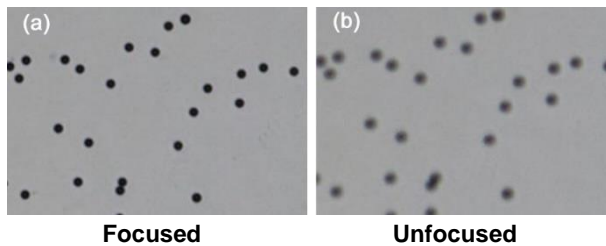
Procedure

1. Invert CCBM-011-2ML vial back and forth 10X
NOTE: DO NOT shake bead solution as it will cause air bubbles to form
2. Vortex CCBM-011-2ML vial for 10 sec on high
3. Invert CCBM-011-2ML vial back and forth 10X
NOTE: DO NOT shake bead solution as it will cause air bubbles to form
4. Remove 20 μL of bead suspension from vial and load into disposable counting chamber



Capillary force automatically spreads cell sample within the counting chamber.

5. Insert slide chamber into Cellometer[®] Automated cell counting instrument
6. Select appropriate Assay and Cell Type Parameters:
 - a. For Cellometer-branded software, assays are included in the database
 - b. For Cellometer K2 using Matrix software, use "Check Bead Assay"
NOTE: If you do not have the appropriate assay, please contact support at Cellc-support@revvity.com
7. Preview Image
8. Adjust focus, as needed



9. Click **Count**
10. Confirm bead counts and viability results are within the lot specific CofA

Store at 2-8°C, protected from light
For Research Use Only

www.revvity.com/cellcountingreagents
Email: Cellc-support@revvity.com