

Cellometer[®] Simply Counted

Reference Beads

Cat #	Bead diameter (µm)	Volume (mL)	Aprox. concentration million/mL
B05-02-050	5	2	5
B10-02-020	10	2	2
B15-02-010	15	2	1

Procedures

1. Invert tube back and forth 10X
2. Vortex tube for 10 sec on high
3. Invert tube back and forth 10X
4. Remove 20µL of bead suspension and add to disposable counting chamber



5. Insert chamber into Cellometer™ automated cell counting instrument



6. Display image
7. Select appropriate Cell Type parameter (see below)

Cat #	Bead diameter (µm)	Cell Type Parameter
B05-02-050	5	Small Cell Size
B10-02-020	10	Medium Cell Size
B15-02-010	15	Large Cell Size



www.nexcelom.com

Tel: 978-327-5340 Fax: 978-327-5341 Email: support@nexcelom.com

Cellometer[®] Simply Counted

8. Click Count
9. Analyze results:
Did the count fall within the expected concentration range?
✓ Yes ... test is complete
✗ NO ... is sample in focus?
✓ Yes ... repeat steps 1 through 9, if still NO then call customer support (978-327-5340).
✗ NO ... adjust focus and repeat steps 1 through 9.

Store at Room Temperature
For Research Use Only



www.nexcelom.com

Tel: 978-327-5340 Fax: 978-327-5341 Email: support@nexcelom.com