

Polystyrene Beads in a Trypan Blue Solution

Materials Supplied

Cat #	Bead diameter (µm)	Total Volume (mL)	Approx. Conc. (x 10 ⁶ beads/mL)
B05-02-050	5	4	5
B10-02-020	10	4	2
B15-02-010	15	4	1

Procedure

1. Invert vial back and forth 10X.
NOTE: DO NOT shake bead solution as it will cause air bubbles to form.
2. Vortex vial for 10 sec on high.
3. Invert 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 a disposable counting chamber.
NOTE: Loading volume for Cellometer™ Ascend™ 8-chamber slide is 10 µL
5. Insert slide chamber into Cellometer automated cell counting instrument.
6. Select appropriate Cell Type parameter:

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

7. Preview image.
8. Adjust focus, as needed.
9. Click **Count**.
10. Confirm bead concentration is within the lot specific CofA.
NOTE: If bead concentration does not fall within the expected range, ensure image is in focus and beads are mostly counted. Contact Revvity Support if assistance is required.

Store at RT (10 to 30 °C)
For Research Use Only

www.revvity.com/cellcountingreagents
Email: USCAN.service@revvity.com