

Counting Cells with a Hemocytometer and a Microscope

1. Prepare a hemocytometer slide. Place a coverslip in the middle of the hemocytometer.
2. In a micro test tube, combine 10 μ l of the cell suspension with 10 μ l of TC10™ trypan blue dye. Gently pipet up and down ten times to mix the cells and dye.
3. Load 10 μ l of the mixture into the opening of either of the two hemocytometer chambers.
4. Expel a droplet of the suspension and let it be drawn under the coverslip by capillary action.

Note: it is important not to overfill or underfill the chamber.

5. Place the hemocytometer in the microscope and allow the cells about 30 sec to settle.
6. Select the 10 \times objective and focus on the grid lines in the chamber. Move the slide so that the field you see is on the central area of the grid (central grid area is 1 \times 1 mm).
7. Count the cells in the four (4) grid areas labeled as 1, 3, 7, and 9. Are the cells in or out? The usual practice is to include cells overlapping the top and left lines but not those overlapping the bottom or right lines – this eliminates redundant counting if adjacent regions are counted.

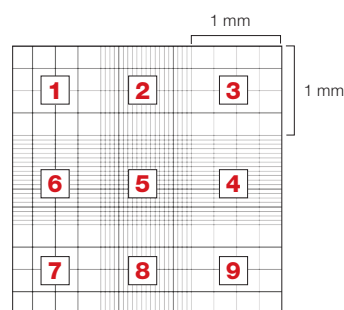


Fig. 1. Neubauer counting grid. The grid is divided in nine 1 mm² sections. Numbers indicate the order in which the sections were counted.

12. Count between 100 and 300 cells
13. Calculate the cell concentration in cells/ml; this is equivalent to the average cell count \times 2,500



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 5044 5699
Canada 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00
France 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 03 963 6050 **Italy** 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666
New Zealand 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04
Singapore 65 6415 3188 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55
Taiwan 886 2 2578 7189 **Thailand** 800 88 22 88 **United Kingdom** 020 8328 2000
