

Biotechnology: A Laboratory Skills Course and Teacher Supplement

Errata and Corrections

Duplication of any part of this document
is permitted for classroom use only.

**IMPORTANT: These corrections are for the
Biotechnology A Laboratory Skills Course
Textbook and Teacher Supplement,
First Edition**

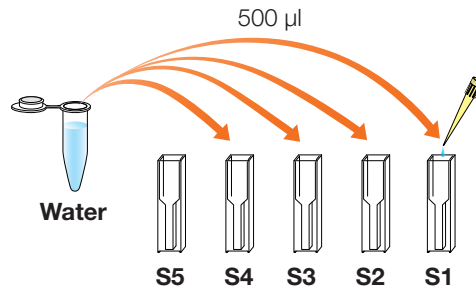
For questions and concerns please contact
Bio-Rad Technical Support at support@bio-rad.com
or by phone at **1-800-424-6723**

Biotechnology: a Laboratory Skills Course, First Edition

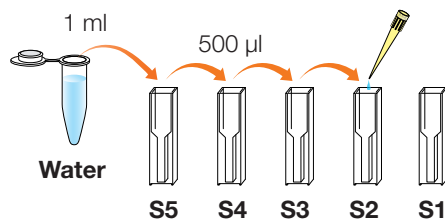
Page 255

Steps 1 and 2 should read:

1. (Optional) Prepare a two-fold serial dilution of *p*-nitrophenol standards. Label five cuvettes **S1–S5**. Add 500 μl of dH_2O to the **S1–S4** cuvettes.



2. (Optional) Add 1,000 μl of 200 nmol/ml *p*-nitrophenol to the cuvette labeled **S5**. Add 500 μl from the **S5** cuvette to the **S4** cuvette and mix. Add 500 μl from the **S4** cuvette to the **S3** cuvette and mix. Add 500 μl from the **S3** cuvette to the **S2** cuvette and mix. Remove 500 μl from the **S2** cuvette and discard. Do not add *p*-nitrophenol to the **S1** cuvette.



Teacher Supplement, First Edition

Page 140

In step 3, the second sentence should read:

Add 100 μl of concentrated enzyme to 25 ml of 1x resuspension buffer and mix by shaking.

Page 141

The Inventory Table should indicate the "quantity per class" of concentrated enzyme as 100 μl .



**Bio-Rad
Laboratories, Inc.**

*Life Science
Group*

Web site bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 1 877 89 01 177 **Belgium** 32 (0)3 710 53 00 **Brazil** 55 11 3065 7550
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 45 44 52 10 00 **Finland** 358 09 804 22 00
France 33 01 47 95 69 65 **Germany** 49 89 31 884 0 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 972 03 963 6050 **Italy** 39 02 216091 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 (0)318 540 666
New Zealand 64 9 415 2280 **Norway** 47 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 (0) 861 246 723 **Spain** 34 91 590 5200 **Sweden** 46 08 555 12700 **Switzerland** 41 026 674 55 05
Taiwan 886 2 2578 7189 **Thailand** 66 662 651 8311 **United Arab Emirates** 971 4 8187300 **United Kingdom** 44 020 8328 2000

