

## Laboratory Quick Guide

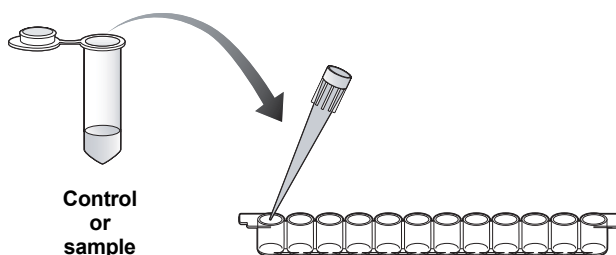
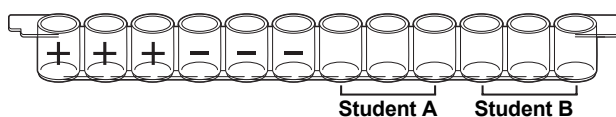
### Antigen Detection ELISA

#### Student Workstation Checklist

One workstation serves 4 students.

Item (Label)	Contents	Number	(✓)
Yellow tubes	Student test samples (0.25 ml)	4	<input type="checkbox"/>
Violet tube (+)	Positive control (0.5 ml)	1	<input type="checkbox"/>
Blue tube (-)	Negative control (0.5 ml)	1	<input type="checkbox"/>
Green tube (PA)	Primary antibody (1.5 ml)	1	<input type="checkbox"/>
Orange tube (SA)	Secondary antibody (1.5 ml)	1	<input type="checkbox"/>
Brown tube (SUB)	Enzyme substrate (1.5 ml)	1	<input type="checkbox"/>
12-well microplate strips		2	<input type="checkbox"/>
50 $\mu$ l fixed-volume micropipet or 20–200 $\mu$ l adjustable micropipet		1	<input type="checkbox"/>
Yellow tips		10–20	<input type="checkbox"/>
Disposable plastic transfer pipet		1	<input type="checkbox"/>
70–80 ml wash buffer in beaker	Phosphate buffered saline with 0.05% Tween 20	1	<input type="checkbox"/>
Large stack of paper towels		2	<input type="checkbox"/>
Black marking pen		1	<input type="checkbox"/>

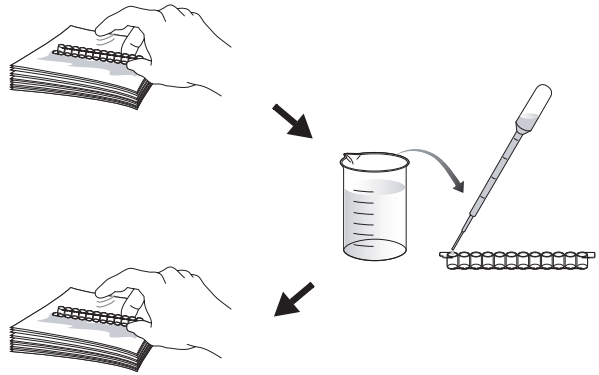
1. Label the yellow tubes with each student's initials.
2. Label your 12-well strip. On each strip label the first 3 wells with a "+" for the positive controls and the next 3 wells with a "-" for the negative controls. Label the remaining wells with your and your partner's initials (3 wells each).
3. Use a fresh pipet tip to transfer 50  $\mu$ l of the positive control (+) into the three "+" wells.
4. Use a fresh pipet tip to transfer 50  $\mu$ l of the negative control (-) into the three "-" wells.
5. Transfer 50  $\mu$ l of each of your team's samples into the appropriately initialed three wells, using a fresh pipet tip for each sample.
6. Wait 5 minutes while the proteins in the samples bind to the plastic wells.



**PROTOCOL II**  
**Antigen Detection ELISA**

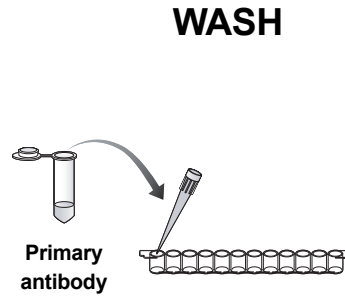
7. WASH:

- a. Tip the microplate strip upside down onto the paper towels, and gently tap the strip a few times upside down. Make sure to avoid splashing samples back into wells.
- b. Discard the top paper towel.
- c. Use your transfer pipet to fill each well with wash buffer, taking care not to spill over into neighboring wells. Note: the same transfer pipet is used for all washing steps.
- d. Tip the microplate strip upside down onto the paper towels and tap.
- e. Discard the top 2–3 paper towels.



8. Repeat wash step 7.

9. Use a fresh pipet tip to transfer 50  $\mu$ l of primary antibody (PA) into all 12 wells of the microplate strip.

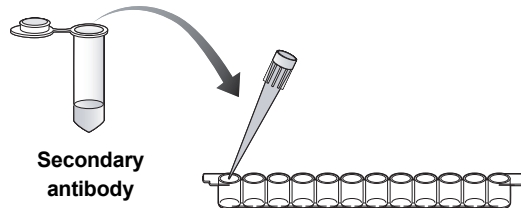


10. Wait 5 minutes for the antibodies to bind to their targets.

11. Wash the unbound primary antibody out of the wells by repeating all of wash step 6 **two** times.

**WASH 2x**

12. Use a fresh pipet tip to transfer 50  $\mu$ l of secondary antibody (SA) into all 12 wells of the microplate strip.

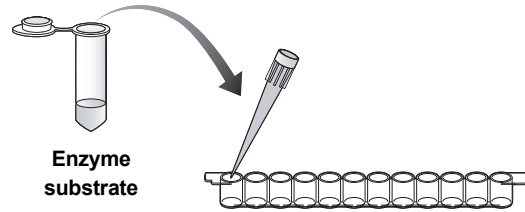


13. Wait 5 minutes for the antibodies to bind to their targets.

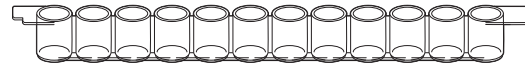
14. Wash the unbound secondary antibody out of the wells by repeating wash step 6 **three** times.

### WASH 3x

15. Use a fresh pipet tip to transfer 50  $\mu$ l of enzyme substrate (SUB) into all 12 wells of the microplate strip.



16. Wait 5 minutes. Observe and record the results.





**Bio-Rad  
Laboratories, Inc.**

*Life Science  
Group*

**Web site** [www.bio-rad.com](http://www.bio-rad.com) **USA** (800) 4BIORAD **Australia** 02 9914 2800 **Austria** (01)-877 89 01 **Belgium** 09-385 55 11 **Brazil** 55 21 2527 3454  
**Canada** (905) 712-2771 **China** (86 21) 6426 0808 **Czech Republic** + 420 2 41 43 05 32 **Denmark** 44 52 10 00 **Finland** 09 804 22 00  
**France** 01 47 95 69 65 **Germany** 089 318 84-0 **Greece** 30 210 777 4396 **Hong Kong** (852) 2789 3300 **Hungary** 36 1 455 8800  
**India** (91-124)-2398112/3/4, 5018111, 6450092/93 **Israel** 03 951 4127 **Italy** 39 02 216091 **Japan** 03-5811-6270 **Korea** 82-2-3473-4460  
**Latin America** 305-894-5950 **Mexico** 55-52-00-05-20 **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** 7 095 721 1404 **Singapore** 65-64153188 **South Africa** 00 27 11 4428508  
**Spain** 34 91 590 52 00 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55 **Taiwan** (886 2) 2578 7189/2578 7241 **United Kingdom** 020 8328 2000