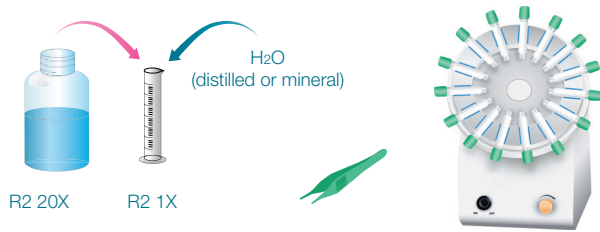


ENOLOGICAL DIAGNOSIS • QUICK GUIDE

# VINEO™ Unstable Proteins 354-8121

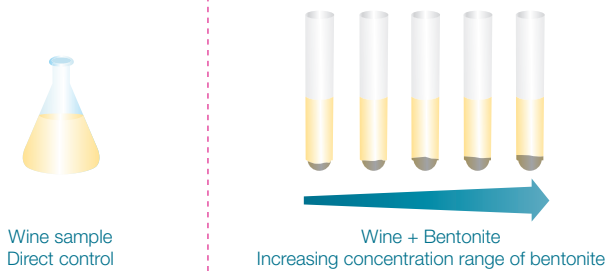


## Equipment Preparation



- Prepare the required number of test strips and incubation chambers
- Dilute the R2 wash solution 1X (40 ml 1X per test strip)
- Hold the solutions at room temperature for 15 minutes before use

## Sample Preparation



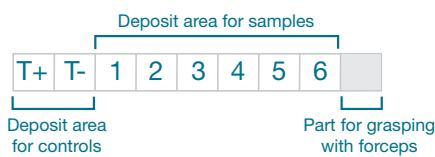
### Direct control of the presence of unstable proteins in wine

- No wine treatment

### Control in wine sample treated with increasing amounts of bentonite

- Prepare a concentration range of bentonite in 2 ml or 4 ml of wine (see annex 1 instructions)
- Mix by inverting 5 to 6 times
- Let precipitate overnight

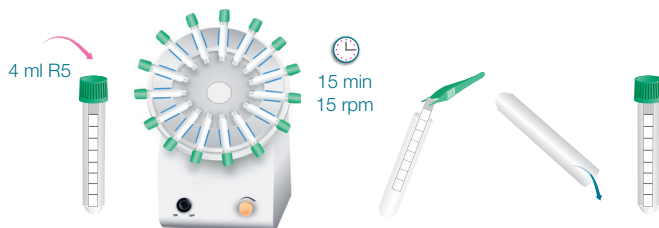
## Sample Deposit



- 5 µl in sample areas 1 to 6

- 5 µl of untreated wine sample in area 1
- 5 µl of the various concentrations within the range of interest into the areas 2 to 6

## Saturation



- Deposit 5 µl of R4 in the area marked T+
- Deposit 5 µl of R3 in the area marked T-
- Let air dry for 10 minutes

- Place the test strip in an incubation chamber
- Add 4 ml of R5 saturation solution
- Place on rotary shaker for 15 minutes at 15 rpm
- Take out the strip using forceps
- Empty the incubation chamber
- Put the strip back into the same tube

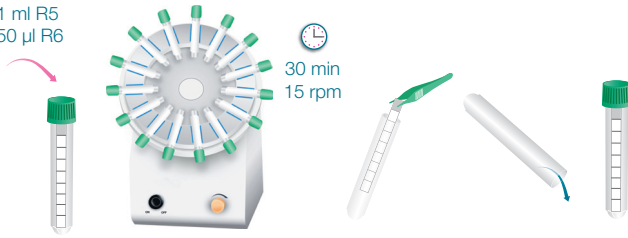
Please read the kit instruction manual for complete and detailed instructions.



## Immuno-detection of Unstable Proteins

### 1 - Primary Detection

3 ml R2  
1 ml R5  
50 µl R6



- Add in the incubation chamber:
  - 3 ml of R2 solution diluted to 1X
  - 1 ml of R5
  - 50 µl of R6
- Place on rotary shaker for 30 minutes at 15 rpm
- Take out the strip using forceps
- Empty the incubation chamber
- Put the strip back into the same tube

### 2 - Washing\*



#### Washing 1:

- Add 4 ml of R2 solution diluted to 1X
- Rinse 10 seconds by manually inverting tube
- Take out the strip using forceps
- Empty the incubation chamber
- Put the strip back into the same tube

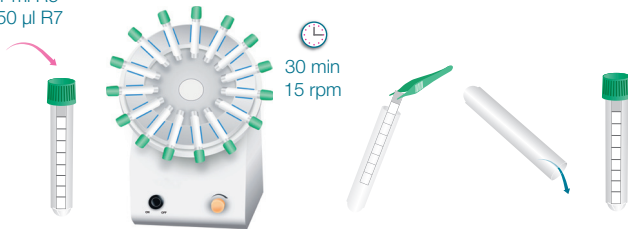


#### Washing 2 (Repeat this step 2 times):

- Add 4 ml of R2 solution diluted to 1X
- Place on rotary shaker for 5 minutes at 15 rpm
- Take out the strip using forceps
- Empty the incubation chamber
- Put the strip back into the same tube

### 3 - Secondary Detection

3 ml R2  
1 ml R5  
50 µl R7



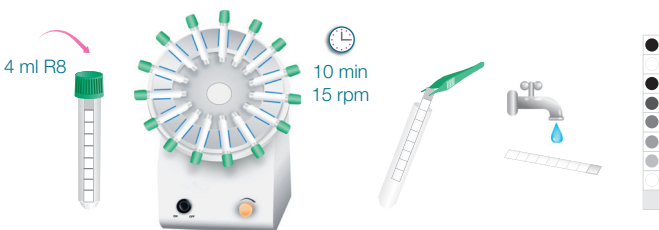
- Add in the incubation chamber:
  - 3 ml of R2 solution diluted to 1X
  - 1 ml of R5
  - 50 µl of R7
- Place on rotary shaker for 30 minutes at 15 rpm
- Take out the strip using forceps
- Empty the incubation chamber
- Put the strip back into the same tube

### 4 - Washing



\* Repeat step: 2 - Washing 1 + 2X Washing 2

## Results



- Add 4 ml of R8 development solution
- Place on rotary shaker for 10 minutes at 15 rpm
- Take out the strip using forceps
- Rinse the strip with water
- Let air dry on a piece of absorbent paper
- Read and interpret results

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://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 31 3689 6600  
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 30 210 777 4396 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 Malaysia 60 3 2117 5260 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 3170 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 66 2 6518311 United Kingdom 020 8328 2000