

VINEO[™] Brettanomy test PCR Kit

Molecular Diagnostics Serving Enology

- Specific: targets the spoilage yeast Brettanomyces bruxellensis
- Safe: controls the risk of ethylphenol production
- Fast: provides results in less than 4 hr
- Quantitative: can be used to quantify the *B. bruxellensis* risk in order to adapt the level of prevention and/or reaction
- Robust: obtains reliable results throughout the process, from fermenting musts through finished wines
- Comprehensive: offers the ability to perform PCR analysis of *B. bruxellensis* and *Zygosaccharomyces bailii* at the same time
- Easy to use: CFX Manager[™] Software IDE automates quantitative and intuitive analysis

Visit us on the web at bio-rad.com/web/VINEO for more information.



VINEO Brettanomytest PCR Kit



VINEO Extract DNA Kit



Detection and Quantification of B. bruxellensis by PCR

B. bruxellensis is a yeast responsible for the growth of 4-ethylphenol and 4-ethylguaiacol, which can impart a strong "off" taste and odor and can affect a wine throughout its development.

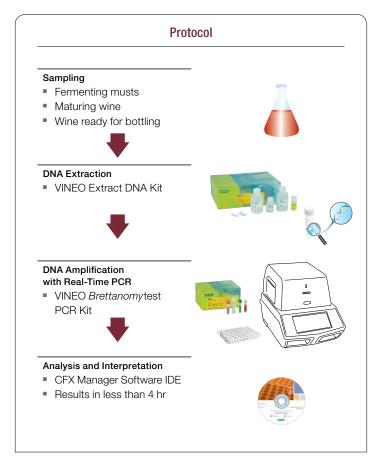
Compared with the traditional microbiological method and other less specific methods, the molecular biology of PCR provides solutions for quick detection with high sensitivity and specificity. VINEO *Brettanomy*test PCR Kit is a quantitative test for the specific detection of *B. bruxellensis* in wines and fermenting musts using real-time PCR.

Implementing this test provides a quantitative result less than 4 hours after DNA extraction using VINEO Extract DNA Kit (catalog #3548100). CFX Manager Software Industrial Diagnostic Edition uses an automatic and quantitative analysis to measure the risk of *B. bruxellensis* in a wine sample. The software then proposes an interpretation of the risk level to the user based on the number of colony forming units (CFU)/ml detected in the wine.

Two DNA extraction protocols are available

Interest	1.8 ml Quantification	45 ml Presence/Absence
Wine in fermentation	•	_
Wine in maturation	•	—
Wine ready for bottling	•	•

The ability to accurately and reliably measure the quantity of *B. bruxellensis* accompanies you in decision-making throughout the winemaking process.



Ordering Information

-	
Catalog #	Description
Kits	
3548100	VINEO Extract DNA Kit, 96 tests
3548101	VINEO Brettanomytest PCR Kit, 96 tests
3548102	VINEO ZYGO bailii Kit, 96 tests
Instruments	
3594990	Industrial Diagnostic CFX96 Touch [™] Deep Well Real-Time
	PCR Detection System, 96 wells (#3600037 for North America)
3593593	CFX Manager Software Industrial Diagnostic Edition
	(#3600042 for North America)
Consumables	
MLL9651XTU	96-Well Low-Profile Unskirted PCR Plates, white, 25
MLL4851XTU	48-Well Low-Profile Unskirted PCR Plates, white, 50
TLS0851XTU	Low-Profile 0.2 ml PCR 8-Tube Strips without Caps,
	thin-walled, white, 120
TCS0803XTU	Optical Flat 8-Cap Strips, for 0.2 ml tubes and plates,
	ultraclear, 120



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 026674 55 05 Taiwan 886 2 2578 7189 Thailand 66 2 651 8311 United Arab Emirates 971 4 8187300 United Kingdom 44 020 8328 2000

(2)