



VINEO™ *ZYGO bailii* PCR Test Kit

Molecular Diagnostics Serving Enology

- Real-time PCR provides high specificity and sensitivity
- Results obtained in less than 4 hr
- CFX Manager™ Software IDE automates qualitative and intuitive analysis
- PCR analysis of *Zygosaccharomyces bailii* and *Brettanomyces bruxellensis* can be done at the same time
- Kits are both easy and ready to use
- Samples extracted using the VINEO Extract DNA Kit can be used to test for both *Z. bailii* and *B. bruxellensis*

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



VINEO ZYGO *bailii* Kit



VINEO Extract DNA Kit



Preventing Re-fermentation of Residual Sugars

Z. bailii is a yeast responsible for significant economic losses in the food industry. Many acidic and/or high sugar content foods are normally considered stable in storage. That is, they readily inactivate a wide range of food-borne microorganisms. However, these same products are susceptible to spoilage by *Z. bailii*.

Its presence is difficult to detect by conventional culture methods, which take a long time and are nonspecific. Rapid, specific, and early detection of this yeast allows producers to take measures to eliminate it before the onset of damage and to prevent unwanted sparkling or the development of any "off" tastes and odors.

Compared with traditional microbiological methods, the molecular biology of real-time PCR provides detection solutions with high levels of speed, sensitivity, and specificity.

Ordering Information

Catalog # Description

Kits

3548100
3548101
3548102

VINEO Extract DNA Kit, 96 tests
VINEO *Brettanomyces* PCR Kit, 96 tests
VINEO ZYGO *bailii* Kit, 96 tests

Instruments

3594990

3593593

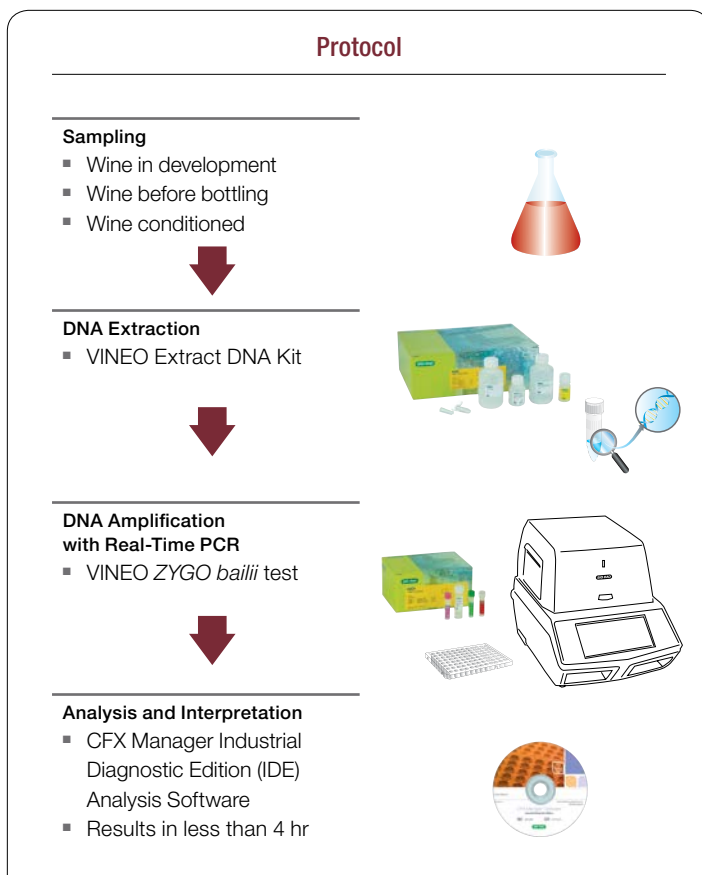
Industrial Diagnostic CFX96 Touch™ Deep Well Real-Time PCR Detection System, 96 wells (#3600037 for North America)
CFX Manager™ Software Industrial Diagnostic Edition (#3600042 for North America)

Consumables

MLL9651XTU
MLL4851XTU
TLS0851XTU

TCS0803XTU

96-Well Low-Profile Unskirted PCR Plates, white, 25
48-Well Low-Profile Unskirted PCR Plates, white, 50
Low-Profile 0.2 ml PCR 8-Tube Strips without Caps, thin-walled, white, 120
Optical Flat 8-Cap Strips, for 0.2 ml tubes and plates, ultraclear, 120



BIO-RAD

**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 2 651 8311 **United Arab Emirates** 971 4 8187300 **United Kingdom** 44 020 8328 2000

