



VINEO™ *ZYGO bailii* PCR Test

PCR Excellence Serving Enology

- Specific: targets only *Zygosaccharomyces bailii*
- Fast: detection in less than 4 hours by real-time PCR
- Sensitive: ensures the absence of “Zygo”
- Easy to use

For more information visit us on the web at
www.foodscience.bio-rad.com



VINEO™ *ZYGO bailii* Test



VINEO™ Extract DNA Kit

BIO-RAD

Prevention of Refermentation of Residual Sugars

Zygosaccharomyces bailii is a yeast responsible for significant economic losses in food industry. Many acids and/or high sugar content such as fruit concentrates, wine, soft drinks, syrups, ketchup, mayonnaise, pickles, salad dressing... are normally considered stable in storage, ie they readily inactivate a wide range of food-borne microorganisms. However, these products are susceptible to be spoiled by *Zygosaccharomyces bailii*. This is a particularly robust species to many stress conditions encountered in food matrices where other microorganisms are not able to grow or even survive. It is described, for its ability to ferment residual sugars in beverages such as wines and fruit juices.

In wines, its presence is difficult to detect by conventional culture methods that are long and non-specific. Rapid, specific and early detection of this yeast in the wine making process can allow to take measures to prevent and eliminate it before the onset of damage which is responsible for olfactory and gustatory deviations, and unwanted sparkling. Faced with traditional microbiological methods, molecular biology provides detection solutions with high levels of speed, sensitivity and specificity.

Key Benefits

- A common DNA extraction for "Brett" and "Zygo": VINEO Extract DNA Kit
- Specificity and sensitivity of real-time PCR
- Result 2 hours after DNA extraction
- Automatic and qualitative analysis: CFX Manager IDE
- Same time PCR analysis possible of « Brett » and « Zygo »
- Easy and ready-to-use kits

Ordering Information

Catalog # Description

VINEO™ Extract DNA Kit
354-8100 96 tests

VINEO™ ZYGO *bailii* Test
354-8102 96 tests

Instruments and consumables

Instruments

359-4990 Industrial Diagnostic CFX96™ Deepwell Real-Time PCR Detection System, 96 wells (360-0037 for North America)
359-3995 Industrial Diagnostic Real-Time PCR Detection System MiniOpticon™, 48 puits
359-3893 CFX Manager™ Software Industrial Diagnostic Edition

Consumables

MLL-9651¹⁻² 96-well low-profile unskirted PCR plates, white, x 25
MLL-4851¹⁻² 48-well low-profile unskirted PCR plates, white, x 50
MSB-1001 Microseal® optically clear adhesive seal, x 100
TLS-0851¹⁻² Low-profile 0.2 ml PCR 8-tube strips without caps, thin-walled, white, x 120
TCS-0803¹⁻² Optical flat 8-cap strips, for 0.2 ml tubes and plates, ultra-clear, x 120



Protocol

Sample

- Wine in development
- Wine before bottling
- Wine conditioned

DNA Extraction

VINEO™ Extract DNA Kit



DNA Amplification with real-time PCR

VINEO™ ZYGO *bailii* Test



Analysis and Interpretation

CFX™ Manager Software IDE

Results in less than 4 hrs



- (1) For items ordered in North America, the following catalog #s apply: MLL-4851XTU, TLS-0851XTU, TCS-0803XTU, MLL-9651XTU
(2) For items ordered in Europe, the following catalog #s apply: 35MLL4851, 35TLS0851, 35TCS0803, 35HSP-9651

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site 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