



TRINEST™ INCUBATOR SHAKER

High quality instrument for high precision results

The 1296-0050 TriNEST™ microplate incubator shaker from PerkinElmer is a high quality device for processing assays on standard-height microplates. Advanced electronic and mechanical design provides the technical excellence you need in order

to obtain quality results in either clinical or research assays.

Compact, versatile solution for plate-based assays

TriNEST can accommodate three plates simultaneously, and with its footprint requires only 388 x 310 mm of laboratory bench space. You can also stack up to three units to give a nine-plate capacity for the same footprint. The instrument is easy to use, with its own keypad and display. Up to 20 incubating and shaking programs can be stored, in which temperature, shaking speed and time are specified. Interval shaking is also supported. You can specify incubation temperatures from 3°C above ambient to 70°C, a very wide temperature range, making the device suitable for most plate-based assays.

Controlled incubation and orbital shaking to assure results quality

To assure highly repeatable results, temperature is accurate to 0.5°C and temperature uniformity across the plate is good, with variation under 0.6°C. For assays requiring simultaneous shaking, the orbital shaker provides an amplitude range from 400 to 1200 rpm, with shaking precision $\pm 2\%$.

Designed for long lifetime and low maintenance

TriNEST has been designed to provide excellent service with materials such as Teflon and acid proof steel to reduce the risk of

corrosion. As well as helping to assure a longer lifetime, careful use of materials maximizes the period between maintenance calls.

SPECIFICATIONS

TriNEST Incubator Shaker

GENERAL

Capacity: 3 plates per unit, SBS standard height
96-well and 384-well plates

Programmable: up to 20 programs

Relative humidity: 15-85 %

LCD graphics display

External power supply

Supports use of plastic or aluminum plate sealers

The 3 plates are processed with the same program but can be loaded at different times

Buzzer and LED lights indicate when incubation is complete

CE marked in conformance with the IVD Directive 98/79/EC

INCUBATION

Temperature range: from 3°C above ambient to 70°C
in 1°C increments

Precision: ± 0.5°C

Variation across plate: < 0.6°C

Warming speed: less than 15 min from +24°C to +70°C

Incubation time: 0-48h in increments of 1 min

Preheat function assuring incubation starts at precisely the specified temperature

SHAKING

Shaking frequency: 400-1200 rpm (or 0 rpm) in
100 rpm increments

Precision: ± 2%

Orbit: ~2 mm

Time: 0-48h in increments of 1 min

Interval shaking: interval up to 60 min in increments of 1s or 1 min

ENVIRONMENTAL CONDITIONS

Operating: 15°C to +35°C, RH 10-75%

Transport: 0°C to + 60°C, RH 10 -80%
packed in transport package

Storage: 0°C to +60°C, RH 10-20%
packed in transport package

PHYSICAL DATA

Weight: ~15 kg

Dimensions: 388 mm W * 310 mm D * 223 mm H

Voltage: 100 – 120 V or 220 – 240 V, 50 / 60 Hz

ORDERING INFORMATION

1296-0050

TRINEST INCUBATOR SHAKER

All PerkinElmer diagnostic products may not be available in all countries. For information on availability please contact your local representative.

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
Phone: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com

PerkinElmer, Inc.
Wallac Oy
PO Box 10
20101 Turku, Finland
Phone: + 358 2 2678 111
Fax: + 358 2 2678 357

ISO 13485
ISO 9001
CMDCAS
ISO 14001
OHSAS 18001



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

©2012 PerkinElmer, Inc. All rights reserved. PerkinElmer is a registered trademark of PerkinElmer, Inc. All other trademarks depicted are the property of their respective holders or owners. PerkinElmer reserves the right to change this document at any time and disclaims liability for editorial, pictorial or typographical errors.