

Instrumentation



for DELFIA® time-resolved fluorescence
and NCS prompt fluorescence assays

Derive maximum benefit

from the DELFIA® technology



- DELFIA® immunoassay kits are available for a wide range of analytes used both in diagnostics and research. Based on time-resolved fluorometry, the DELFIA method is today's leading non-radioisotopic alternative to RIA.
- To help you to get the best possible results from your DELFIA assays, PerkinElmer provides a complete, optimized instrumentation system for semi-automatic performance of time-resolved fluorometry assays.



A complete optimized system

For reagent dispensing

- The DELFIA® Plate Dispense and DELFIA Dispense Unit (DELFIA assays also work well with manual pipetting as well as with automatic pipetting stations)

For incubation

- The DELFIA Plateshake
- NCS Incubator

For washes

- The DELFIA Platewash

For assays with blood spot samples

- The Wallac DBS Puncher + MultiPuncher
- You can also combine two stages of blood spot assays with the new time-saving DELFIA Washer-Diskremove

For measurement

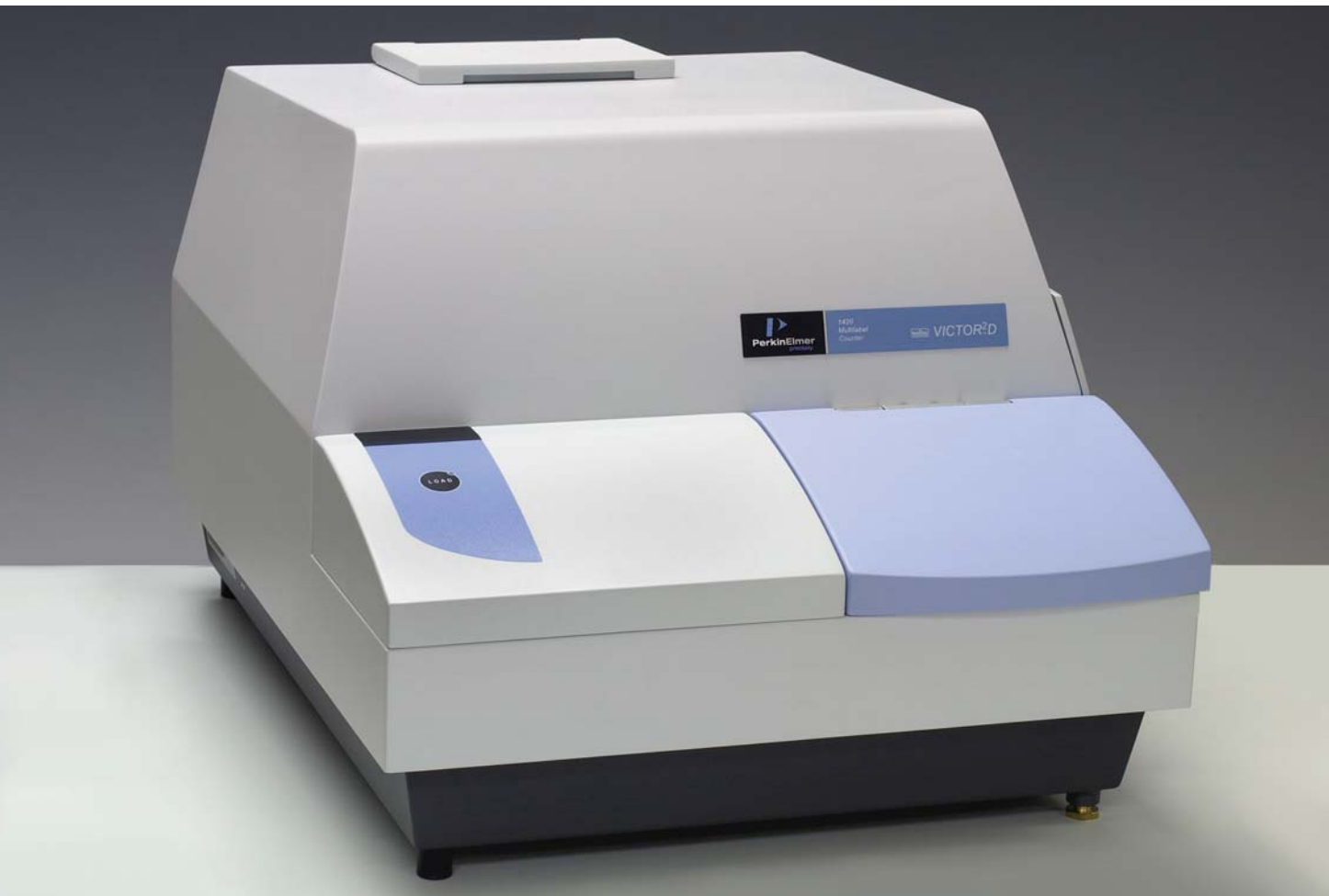
- VICTOR²™ D Fluorometer (or for research use, the VICTOR³™ multilabel counter)

For quality control and data management

- MultiCalc® software

For information about fully automatic performance of DELFIA assays, please ask for details of the Wallac AutoDELFLIA® automatic immunoassay system.

For all DELFIA® assays



Wallac VICTOR²™ D Fluorometer or VICTOR³ multilabel counter

The Wallac VICTOR² D fluorometer is designed for clinical use and is supplied with the ready protocols for all PerkinElmer diagnostic and screening assays based on either time-resolved fluorescence or prompt fluorescence. VICTOR² D includes MultiCalc software for assay evaluation and quality control.

The Wallac VICTOR³ multilabel counter is suitable for a wide variety of research applications. It measures time-resolved fluorescence from four different lanthanides. It is also available in versions that measure assays based on prompt fluorescence, fluorescence polarization, luminescence and absorbance.

DELFIA® Plate Dispense and DELFIA® Dispense Unit

The 1296-041 DELFIA® Plate Dispense allows the automatic and precise addition of Enhancement Solution to all wells or just selected strips. Materials and speeds are optimized for use with DELFIA reagents, to minimize the background and avoid foaming.

Used either alone or with the DELFIA Dispense Unit, the Plate Dispense greatly reduces assay “hands-on” time. The microtitration plate is simply loaded onto the plate holder and the pump is activated. When the plate is ready, the next one can be loaded.

The Plate Dispense can be fitted with the 1296-043 DELFIA Dispense Unit which allows automatic dispensing of buffer or tracer solution. The dispense unit has a completely separate tubing system to avoid contamination. A special rinse function makes it easy to change the liquid being used.

DELFIA® Plateshake

The 1296-003/004 DELFIA Plateshake is a general purpose microtitration plate shaker especially optimized for use with DELFIA diagnostic kits. As well as two set shaking speeds which have been optimized to give the best results for DELFIA assays, you can also select any other speed within the range 100 - 1350 rpm. The constancy of the speeds is controlled by means of an electronic feedback circuit. A rubber mat keeps the plates firmly in place and ensures easy loading.

DELFIA® Platewash

The 1296-026 DELFIA Platewash is a general purpose microtitration plate and strip washer especially standardized for use with DELFIA diagnostic kits.

Control is by softkeys and the operations available are displayed on a two-row, clear, 40-character liquid crystal display. Protocols for all DELFIA kits are provided with the instrument. There is alternatively memory space for up to 75 freely programmable protocols. Dispensing and aspiration probes are paired to avoid overfilling of wells.

The system is self-contained with its own vacuum pump. For even dispensing, a membrane pump is built in. Also supplied are three bottles for wash solution, distilled water and waste. To keep the instrument in a permanent state of readiness, there is an automatic rinse function, and for easy servicing, complete diagnostic software is included.

NCS Incubator

The 1296-008 NCS Incubator has a three-plate capacity and an extended temperature range up to nearly 70°C. Temperature variation is less than 0.3° C across the microplate. To eliminate temperature gradients and edge effects, microplates are evenly heated from all sides. This ensures assay reproducibility.



DELFIA® Plate Dispense and DELFIA® Dispense Unit



DELFIA® Plateshake



DELFIA® Platewash



NCS Incubator

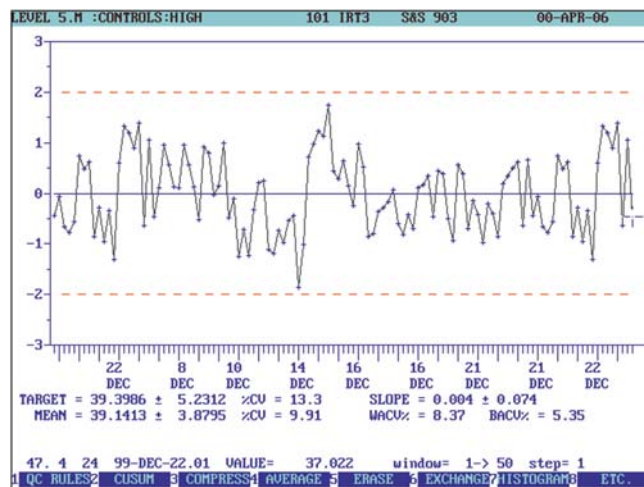
MultiCalc® software

MultiCalc software provides unbeatable data evaluation and quality control for DELFIA® and other biospecific label assays. With MultiCalc, you can handle up to 6 different levels of controls in each assay.

Controls can be stored from 150 different assays to form a Levy-Jennings graph. This gives the user the facility to create a reliable and automatic QC follow-up of assays. Precision profiles and histograms complete the QC part of the program.

With MultiCalc's unique programmable output feature, result reporting can be tailored to the customer's needs. For example, when working with two different assays as part of the same investigation, the results can be combined to provide a clinically meaningful report.

MultiCalc can be connected to a laboratory information system for patient file management. Patient records can then be drawn from the mainframe, updated with the results of DELFIA tests, and then returned to the mainframe.



For DELFIA® bloodspot assays



DELFLA® Washer-Diskremove



Wallac MultiPuncher

DELFLA® Washer-Diskremove

The 1296-0010 DELFLA® Washer-Diskremove is optimized for use with DELFLA neonatal screening assays run on filter paper. It automatically removes the eluted filter paper disks from the wells of the microplate, and also performs all needed wash stages as specified in the assay protocol. Up to 100 protocols may be stored.

The instrument consists of two units, the Washer-Diskremove unit and the separate vacuum unit with a waste bottle.

Wallac MultiPuncher

The 1296-081 Wallac MultiPuncher is an advanced and convenient instrument for punching dried blood spots and dropping them into microplates. Capacity for six plates allows handling of the highest workloads with minimum hands-on time. There are two different sized punch heads (3.2 and 6 mm), and both can be used in the same run. Windows-based software controls punching making for easy operation with minimal user intervention. Wallac MultiPuncher features a bar code reader to ensure positive sample identification and maximum security. A light guide facilitates positioning of the blood spot for optimal punching and the system automatically verifies that the disk falls into the well.



Wallac DBS Puncher

The 1296-071 Wallac DBS Puncher automatically punches dry bloodspot samples into microtitration plates. Three different sizes of bloodspot may be punched (3.2, 4.7 or 6 mm).

Two plates may be loaded simultaneously and the plate height allowed is adjustable, so deep well plates can be accommodated. After punching, disk detection ensures that the disk is correctly placed in the microtitration well.

The DBS Puncher may be connected to a barcode reader for identification of the patient sample. It may be run from an external computer, receiving a plate layout which specifies the sample codes for the wells and the punch sequence.

The instrument is controlled by hand or by footswitch. Ease of use is enhanced by a clear double line display, and by illumination of the area punched.

A complete set of optimized products

1420-020 VICTOR²™ D Fluorometer

Measurement mode	TR-fluorescence, prompt fluorescence
Measurement time	1 second per sample, 3 minutes per plate
Display and peripheral equipment	Includes MultiCalc for assay evaluation and quality control
Barcode reader	Ready protocols for all Wallac analytes Optional barcode reader for worklist identification
Physical dimensions	Height 380 mm (510 mm with stackers), width 485 mm, depth 590 mm, weight 45 kg (1420-020), 56 kg (1420-021 with stackers)

1296-041 DELFIA® Plate Dispense

Speed	3 min 20 sec for one plate
Volume dispensed	200 µl per well (preadjusted)
Accuracy	Better than 1 % at 200 µl
Precision	Better than 3 % at 200 µl
Dimensions	Height 268 mm, width 290 mm, depth 300 mm, weight 10.3 kg.

1296-043 DELFIA® Dispense Unit

Speed	2 min 20 sec for one plate
Volume dispensed	50 or 100 µl per stroke (user adjustable)
Accuracy	Better than 1 % at 200 µl
Precision	Better than 3 % at 200 µl
Dimensions	Height 85 mm, width 150 mm (+ 65 mm for the bottle holder), depth 135 mm, weight 2 kg

1296-003/004 DELFIA® Plateshake

Speed	1 variable setting within the range 100-1350 rpm
Capacity	4 microtitration plates
Shaking amplitude	3 mm
Drive	Split-pole motor
Dimensions	Height 128 mm, width 245 mm, depth 285 mm, weight 5 kg.

1296-026 DELFIA® Platewash

Number of wash methods	10
Number of strips	1-8
Dispensing/aspirating manifold	8, 2 * 8, 3 * 8, 12 or 2 * 12 channels
Wash programs	Special program for each DELFIA kit Alternatively 75 freely programmable
Dispensing	By built-in pump with aspiration to prevent overfilling
Aspiration	By built-in pump
Bottles	3 x 2 litre plastic bottles
Rinse	Automatic rinse function
Voltage / Power	110/220 V 50/60 Hz / 100 W
Dimensions	Height 193 mm, width 325 mm, depth 525.5 mm (plate carrier out), weight 9.6 kg.

1296-008 NCS Incubator

Temperature	up to 70° C
Capacity	3 plates
Dimensions	Width: 400 mm, height: 450 mm, depth: 200 mm

1296-0010 DELFIA® Washer-Diskremove

User interface	2 x 20 characters LCD screen (backlit), 5 diaphragm keys on a flat keypad
Vacuum	Works with an external vacuum unit
Number of wash methods	10
Number of strips	1-8
Manifold	8 or 12 channels
Wash programs	Special program for each DELFIA kit plus 75 freely programmable to total of 100
Bottles	Waste: 8 L, Wash: 5 L, Rinse: 2L
Dimensions	Width: 368 mm, Length: 530 mm (plate carrier out), Height: 483 mm, Weight: 16 kg
Vacuum unit	Width: 140 mm, length: 400 mm, height: 370 mm, weight: 13 kg (approx.)

1296-071 Wallac DBS Puncher

Punch diameter	3.2 mm, 4.7 mm or 6 mm
Plate height	15 - 41 mm
Operational control	2-row, 24 character per row display, 4 selection buttons
Dimensions	Height 305 mm, width 420 mm, depth 480 mm, weight 20 kg

1296-081 Wallac MultiPuncher™

Punch diameter	3.2 mm (1/8") and/or 6 mm
Plate height	15 mm
Input/Output connections	Instrument is connected to the PC via RS-232 Barcode reader is connected to the PC via RS-323
Dimensions	Height 55 cm, width 100 cm, depth 55 cm, weight 45 kg

1420 VICTOR³™ Multilabel counter

Light sources	Continuous light source for fluorometric and photometric measurements Optional flash light source for TR-fluorometric measurements
Detection units	Photometry, Fluorometry, luminometry, TR-fluorometry and fluorescence polarization: depending on the model
Physical dimensions	Height 380 mm (510 mm with stackers), width 485 mm, depth 590 mm, weight 45 kg (1420-012), 56 kg (1420-015 with stackers)

All PerkinElmer diagnostic products may not be available in all countries.

For information on availability please contact your local representative.

PerkinElmer Life and Analytical Sciences

710 Bridgeport Avenue
Shelton, CT 06484-4794 USA
Phone: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com

Wallac Oy, PO Box 10
20101 Turku, Finland
Phone: + 358 2 2678 111
Fax: + 358 2 2678 357
www.perkinelmer.com



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

©2005 PerkinElmer, Inc. All rights reserved. DELFIA and PerkinElmer are registered trademarks of PerkinElmer, Inc. All 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.