

HUMAN HEALTH

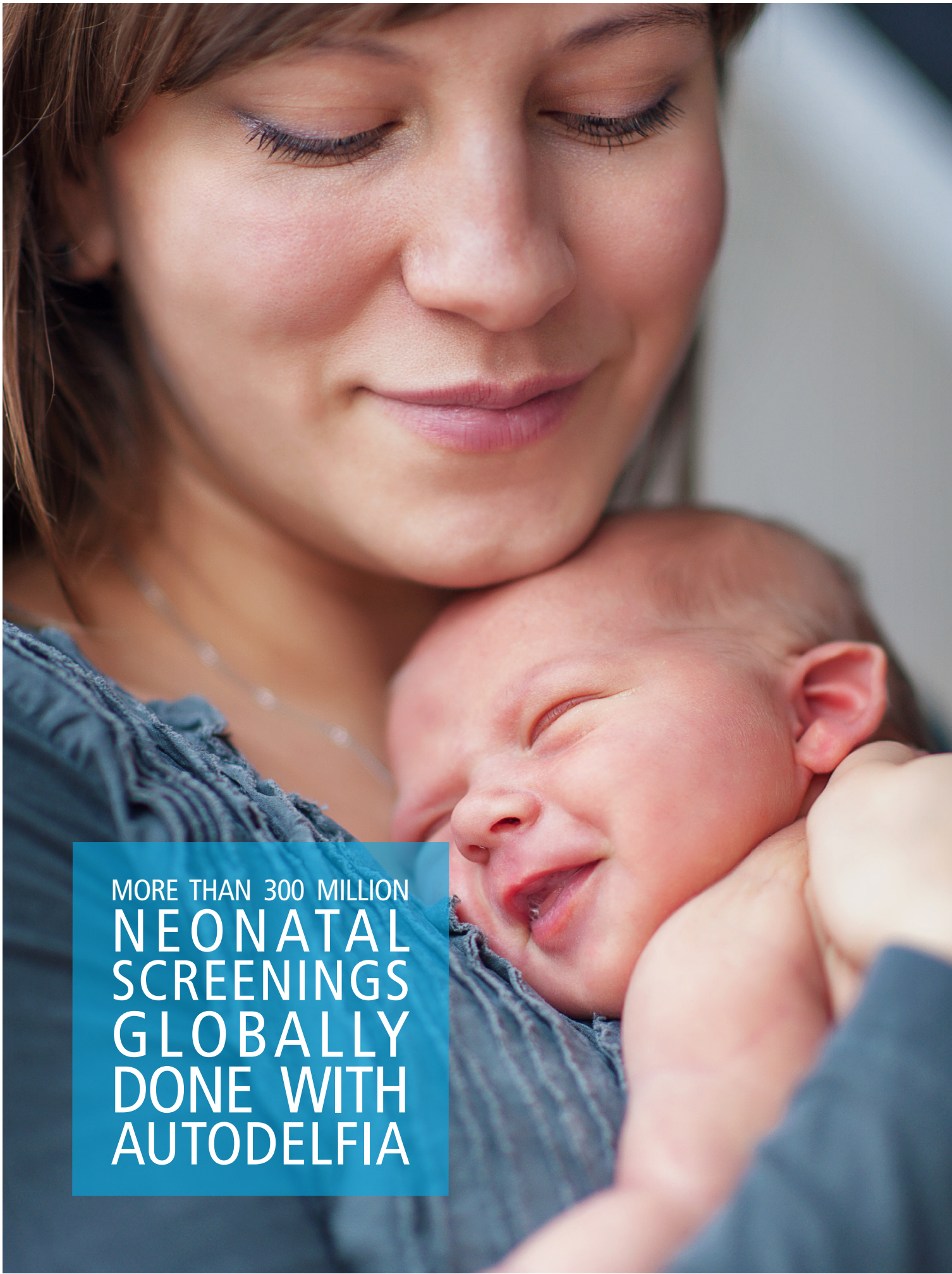
ENVIRONMENTAL HEALTH

MAKE THE CHANGE
FOR BETTER
NEONATAL
SCREENING



Introducing the new AutoDELFIA® immunoassay system
for newborn screening


PerkinElmer
For the Better



MORE THAN 300 MILLION
NEONATAL
SCREENINGS
GLOBALLY
DONE WITH
AUTODELFIA



THE WORLD'S LEADING SCREENING PLATFORM

PerkinElmer's AutoDELFIA immunoassay system is the leading neonatal screening platform with more than 300 million screenings to its credit. Fully automatic and easy to use, AutoDELFIA has become the screening solution of choice for laboratories around the world, and has helped thousands of children start life with a fighting chance.

Now PerkinElmer is proud to introduce the new AutoDELFIA system. By making AutoDELFIA even more effective and easier to use, we give you a clear reason to change your screening for the better.

WHY CHOOSE THE NEW AUTODELFIA

- ✓ Better design for improved user-friendliness
- ✓ Better software for more efficient workflow
- ✓ Better efficiency for enhanced cost-effectiveness.



THE TRUSTED SYSTEM FROM INSIDE TO OUTSIDE

Proven screening record

AutoDELFI is a fully automatic system for sensitive and reproducible quantitation of analytes commonly used in neonatal screening programs.

All kits employ the sensitive and robust DELFIA® chemistry, based on time-resolved fluorescence. Every assay stage including measurement is performed automatically and results are easy to interpret and manage.

Worklists from the punching device can be interpreted automatically, and positive sample IDs are provided for superior traceability if using barcoded sample collections cards.



Optional Sample Processor

AutoDELFI can be equipped with an advanced Sample Processor for serum based screening.

MARKEDLY IMPROVED EFFICIENCY

Designed to bring a new level of convenience and efficiency to neonatal screening.



User-friendly design

Improved ergonomics ensure better working conditions for laboratory personnel.



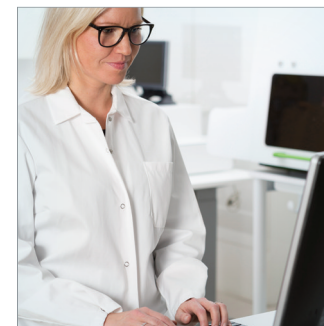
Safer screening

New enclosed plate loading minimizes the risk of contamination.



Laboratory friendly

The front lid moves up instead of out for more convenient instrument access. LED lights inside improve visibility.



Secure and reliable audit trail

Instrument is protected via safe user identification. All change actions are logged to the audit trail for reliable follow-up.



Clean and fast

Color-coded bulk liquid bottles with new super-quick connectors are easy to change and fill up.



Trusted AutoDELFI kits

The AutoDELFI kit includes reagents, QC material and specific QC certificates. Barcoded plates confirm that plates are loaded in correct order, and worklists are picked up based on the plate barcode.



Hands-on convenience

The new instrument table and optional trolley for bulk liquid bottles make for easier handling.



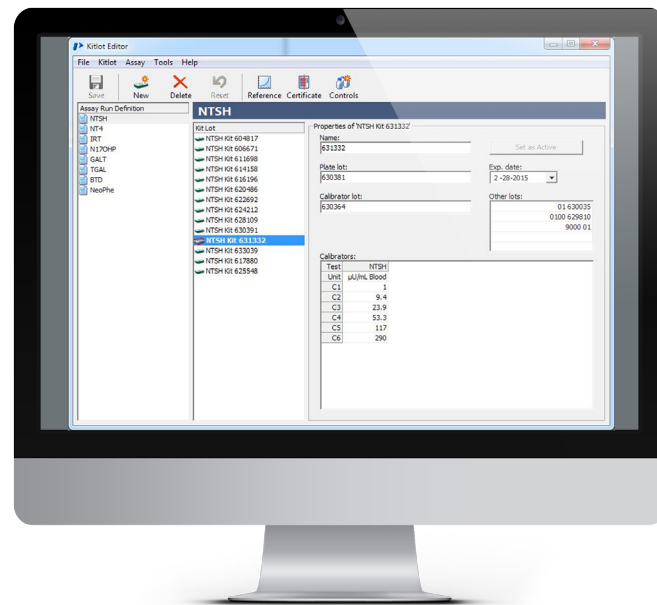
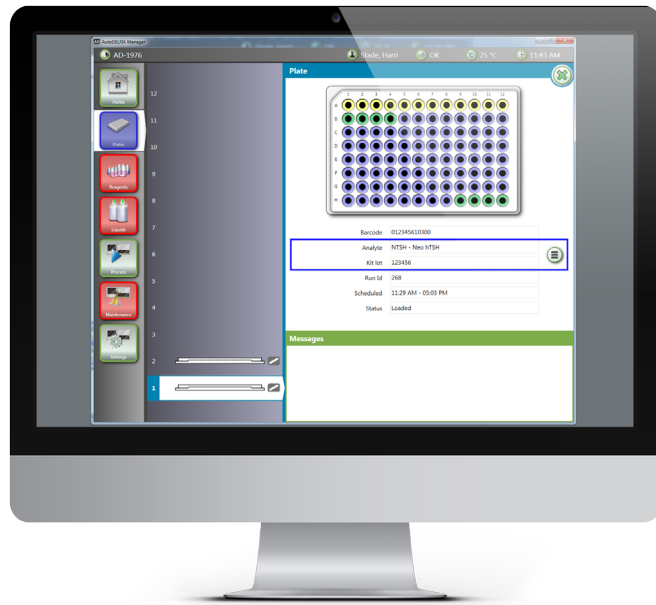
Reliably cost-effective

Robust build and excellent adaptability enhance cost-effectiveness throughout the screening process. PerkinElmer instrumentation and reagents are manufactured to the highest quality and performance standards.



For a more detailed description, please check out the AutoDELFI video at <https://perkinelmer.box.com/AutoDELFI-video>

SMARTER SOFTWARE



AutoDELFI features a completely new Manager instrument user interface program, together with a Workstation program for managing and reporting results

AutoDELFI MANAGER 3.0

Performs essential device operations.

- New graphical user interface and operational modes for improved workflow management
- Easier to learn and use – reducing the need for training
- Fully integrated with the PerkinElmer punchers – worklists are automatically picked up when plates are loaded
- One-glance overview of screening process with new HOME PAGE feature
- Tracking of mandatory maintenance routines by new SYSTEM MAINTENANCE feature
- More effective tracking and management of plates with new PLATE VIEW feature
- Software protection with user specific access rights and audit trail
- Improved reagent pipette error handling and recovery features

AutoDELFI WORKSTATION 3.0

Manages and reports screening results.

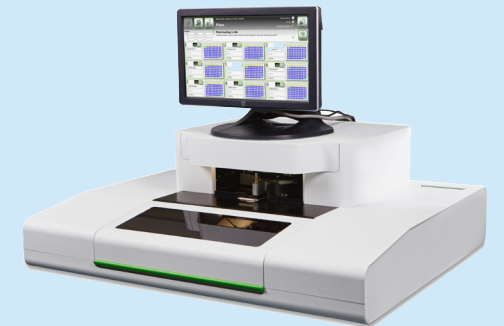
- Replaces the current counts-to-concentration software
- User interface and usability based on SpecimenGate® – PerkinElmer's dedicated informatics solution for newborn screening laboratories
- Includes the same four modules as in the GSP® Workstation: the Kitlot Editor, the Plate Generator, the Result Viewer, and the Quality Control Program
- Automatic import of kit lot data using new 2D-barcode AutoDELFI kit QC certificates
- Improved safety and reliability – no manual work stages and minimal risk of user error



IN GOOD COMPANY

The AutoDELFI system is easy to interface with other PerkinElmer screening devices and software for enhanced screening performance.

1



PUNCHER

Panthera-Puncher™ 9 is a next-generation dried sample punching device that can prepare up to 9 microtitration plates simultaneously. The punch pattern is automatically adapted to the shape and size of the blood spot for more efficient use of sample material without requiring the user to adjust the pattern to fit the area to be punched.

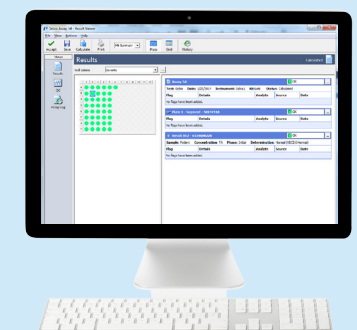
2



PLATE PROCESSOR AND REAGENT KITS

The AutoDELFI Plate Processor is an immunoassay system widely used in newborn screening laboratories. The new AutoDELFI uses the same proven kits as before for screening of some of the most common neonatal screening disorders. The processor performs all assay stages and measurement automatically.

3



SPECIMEN GATE®

Specimen Gate® modules provide control and monitoring tools for the entire screening process from overall screening management to electronic reporting of results and follow-up of screen-positive patients.

HEALTHIER FUTURE

NBS disorders screened by the AutoDELFI System are

- ✓ Congenital Hypothyroidism
- ✓ Congenital Adrenal Hyperplasia
- ✓ Cystic Fibrosis

INSTRUMENT SPECIFICATIONS

Physical dimensions (incl table but excl PC)

Height	1740 mm (2010 mm with lid open)
Width	1246 mm
Depth	910 mm
Weight	270 kg (216 kg without table)

Power requirements

Power consumption max 600VA
Mains voltage 100 - 240 V, 50 - 60 Hz

Environmental conditions

Operating temperature 15 - 30 °C
Relative humidity 20 - 80 %
Noise < 66 dB

Water requirements for wash and rinse

CLRW/CLSI or equivalent deionized water

Samples

Blood disks from dried blood spots punched into 96-well plates (Liquid sample tubes if using the optional Sample Processor)

Loading

Plates, reagents and bulk liquids loaded in a batch
One additional loading of plates, reagents and bulk liquids possible during run

PERFORMANCE SPECIFICATIONS

Capacity

Sample capacity 12 plates (1152 wells)
Reagent capacity up to 4 reagent cassettes (neonatal assays)
Manual filling of wash solution and rinse, no intervention needed during run
Automatic emptying of waste bottle (optional)

Precision

CV% of one plate with 1 nM Eu solution < 1%

Measurement unit

Measurement mode TRF (filters for Eu and Sm)
Linearity error < ±10% up to 6 nM Eu solution (approx. 6 million counts)

Temperature control

Inside temperature controlled to +25 ± 3 °C

ORDERING INFORMATION

1235-5110 AutoDELFIA Plate Processor system

The AutoDELFIA Plate Processor system includes the AutoDELFIA Plate Processor and an external PC with Windows, the AutoDELFIA Manager user interface software, and the AutoDELFIA Workstation software, including the Kitlot Editor, Plate Generator, Result Viewer, and Quality Control modules, with user manuals. A table for the Plate Processor and bottles for bulk liquids are also included.

The AutoDELFIA Plate Processor consists of all modules needed to perform an automatic AutoDELFIA DBS assay: Barcode reader for plates and reagent cassettes, Reagent dispenser, Shaker/Incubator, Disk remover, Washer, Enhancement Solution dispenser, Time-resolved fluorometer, and Plate transport system.

The AutoDELFIA Plate Processor system includes:

- 1235-0110 AutoDELFIA Plate Processor
- 1235-4230 AutoDELFIA Plate Processor table
- 1235-8100 AutoDELFIA V3.0 PC, including the Wallac Instrument PC with LED touch screen, PCI Arcnet card, Windows 7 (or newer), and the 1235-8090 AutoDELFIA V3.0 Software package

Options (sold separately)

- AAAJA-0002 APC Smart-UPS 2200 VA, Uninterruptible Power Supply (230 V)
- UPSAP200 APC Smart-UPS 2200 VA, Uninterruptible Power Supply (115 V)
- 2011-0030 Wallac laser printer (230 V)
- 2011-0040 Wallac laser printer (115 V)
- 1235-4140 Waste pump option
- 1235-4240 Bottle trolley

Kits

- B005-204 AutoDELFIA Neonatal IRT kit (4 plate)
- B005-212 AutoDELFIA Neonatal IRT kit (12 plate)
- B015-112 AutoDELFIA Neonatal 17-OHP kit (12 plate)
- B024-104 AutoDELFIA Neonatal 17α-OHP kit (4 plate)
- B024-112 AutoDELFIA Neonatal 17α-OHP kit (12 plate)
- B032-312 AutoDELFIA Neonatal hTSH kit (12 plate)
- B065-112 AutoDELFIA Neonatal Thyroxine (T4) kit (12 plate)

Consumables (sold separately)

- 1235-411 Dilution vessels (100 pcs)
- 1235-402 Reagent tips (960 pcs)
- B117-100 DELFIA/AutoDELFIA Wash Concentrate (8 x 250 mL)
- B118-100 DELFIA/AutoDELFIA Enhancement Solution (8 x 250 mL)

Products in this brochure may not be available in all countries. Please check availability with your local PerkinElmer representative. The AutoDELFIA Plate Processor system is for use in newborn screening.

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (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

Copyright ©2015, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

1599-9800-01 BRO AutoDELFIA Plate Processor, October 2015 PKI