

HUMAN HEALTH

ENVIRONMENTAL HEALTH

LOOK INTO THE
FUTURE!



GSP

2021 Genetic Screening Processor

Brochure not for distribution in the USA.


PerkinElmer®
For the Better



THE PERFORMANCE
YOU NEED

GSP™ addresses the need for more efficient and more reliable neonatal screening processes. To secure the performance of your operation now and in the future, the new instrument provides greater flexibility and ease of use, reduced hands-on stages and reduced risk of error.

From plate loading to results output, GSP performs all assay stages automatically and at the optimum time. It is used with PerkinElmer dried blood spot assays including those employing the high sensitivity DELFIA® chemistry. It also suits simpler fluorescence or absorbance-based measurement technologies.

For laboratories wishing to implement full automation from specimen reception to results delivery, the GSP instrument is compatible with PerkinElmer's AutoPuncher™ sample processor and Specimen Gate® informatics.



GSP is operated via a built-in touch-screen; actions relating to workflow, including results reception, are performed at a separate PC.



WHAT GSP WILL MEAN IN YOUR LABORATORY

High performance automation – for less hands on time

The GSP system provides automation of every assay stage so that laboratory personnel have more time for other important tasks.

Continuous loading of samples – for flexible workflows

The GSP instrument provides the convenience of continuous sample loading. The operator simply replaces the existing magazine with one containing the new plate or plates. GSP will automatically uptake the loaded plates and schedule the assays.

Barcoding of all materials – for reduced risk of error

To support improved traceability and ease of use, all plates, and all reagents, QC materials and lot specific QC certificate information are barcoded.

Direct access water and waste lines – for further work savings

GSP can be connected directly to a non-ionized water supply and waste-line, so you no longer need to add water and drain the waste

Superior multi-technology solution – for results confidence

PerkinElmer neonatal assays employing prompt fluorescence can now be automated for running on the GSP system.

GSP versions of DELFIA assays utilize a new chelate chemistry with better tolerance towards interfering substances such as EDTA. In addition, for the assay's enhancement stage, DELFIA Inducer assures a superior signal.

GSP™ Automatic System



Reagent storage compartments are temperature controlled so reagents can remain safely in the instrument after the particular assay is complete.



Using the touch screen of the GSP User Interface you can perform the basic operations like adding reagents and emptying waste.

Easy to use – supporting staff rotation

GSP is easy for laboratory staff to learn to use. The instrument is controlled from the touch-screen of the GSP User Interface. This prompts the operator whenever it is necessary to perform an action such as adding reagents or removing plates.

Purpose-designed informatics – for operational efficiency

PerkinElmer informatics solutions are the result of years of experience as the world leader in data management solutions for screening. For use with GSP, new GSP Workstation software is designed to follow laboratory workflow to obtain reliable, quality controlled results with maximum efficiency.

Integrated screening operation – supporting growth and change

Fully compatible with the GSP Workstation product, PerkinElmer's Specimen Gate LIMS provides a backbone for integrated operation of laboratory instruments. The GSP instrument can thus be linked to the AutoPuncher instrument or other punching devices.

Use together with AutoPuncher allows a laboratory to extend automation to include all sample preparation stages from the reception of the dried blood spot cassettes at the laboratory.



SEE HOW EASY YOUR WORK BECOMES

GSP supports operator
comfort and ease-of-use



1 Log in

You operate the instrument from the GSP User Interface, through which you also access all the main functions. As a recognized user, your work with the system will start when you log on and supply your PIN code.

2 See the information you need

The touch screen allows you to see instantly the status of consumables, reagents, and waste capacity.



3 Obtain more detailed information

Press any one of the areas to see the status of the marked item.



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 Genetic Screening
Center of Excellence**
Wallac Oy, PO Box 10
20101 Turku, Finland
Phone: + 358 2 2678 111
Fax: + 358 2 2678 357



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

©2009 PerkinElmer, Inc. All rights reserved. GSP and Wallac are trademarks and 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.