

Punching for the next generation



Introducing the 1296-091 Wallac AutoPuncher™



The first fully automated dried sample punch for the newborn



The Wallac AutoPuncher™ has been designed and developed to meet the growing need for automated tools for dried sample handling in neonatal screening laboratories.

The world's first fully automated dried sample punch instrument, it links to Specimen Gate® and LifeCycle™ LIMS programs to create a fully integrated system.



Wallac AutoPuncher™ frees up operator time by performing all of the routine tasks automatically

The AutoPuncher user simply loads microplates and magazines containing specimen cassettes into the instrument and sets up the punching run with the Workstation software. The Workstation software is easy to use, with every task being clearly explained.

The instrument then performs all operations automatically, normally alerting the user only when it is time to collect the ready prepared microplates and unload the magazines.

Alert module attracts the operator's attention if an action is required

Whether just one AutoPuncher or several units are running, there is no need for watching and waiting. Operators can freely turn their attention to other tasks confident that the system's Alert module will attract their attention should an operator activity be needed.

screening laboratory



Takes 300 specimen cassettes

Wallac AutoPuncher™ accepts 300 specimen cassettes at a time. The instrument checks the ID of the cassette, communicates bi-directionally with the LIMS software, and determines what tests have to be processed and how many spots need to be punched.

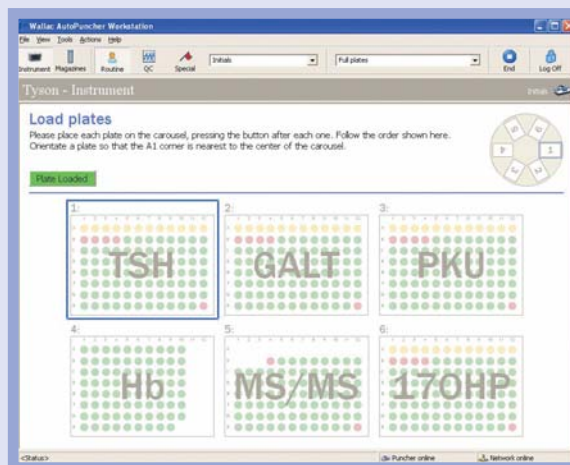
A scanning device automatically checks that there are enough specimens present and sets the punching coordinates. The specimen cassette is then moved to the punching head and the appropriate numbers of samples are punched into their respective plates/trays. The cassette is then returned to the same holding position for either sample follow up or specimen banking.

Checks the condition of the blood spots

As part of its routine operation AutoPuncher scans the dried blood specimen areas on specimen cassettes. It thus ensures that the specimens are suitable for use and determines optimum positions for punching.

Will support robot integration

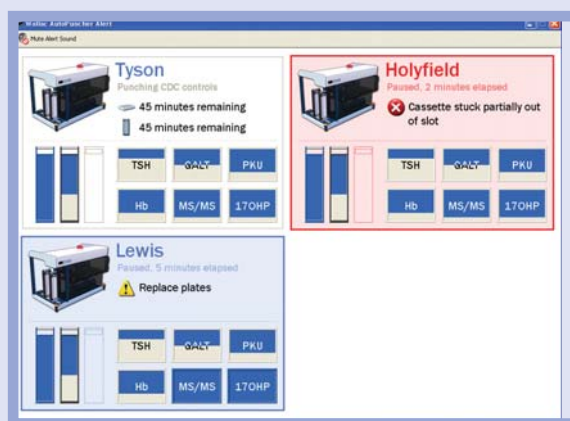
There are six plate loading bays, and each will accommodate one standard microtitration plate/tray or deep well plate. The unique plate holder has been specifically designed for future integration with external plate handling robots. It allows easy loading and unloading of the plate or tray by external plate handling robots.



Running the Wallac AutoPuncher is simple, with the Workstation software guiding the operator at every step. Routine operations require only minimal interaction, allowing more experienced operators to work quickly and efficiently.

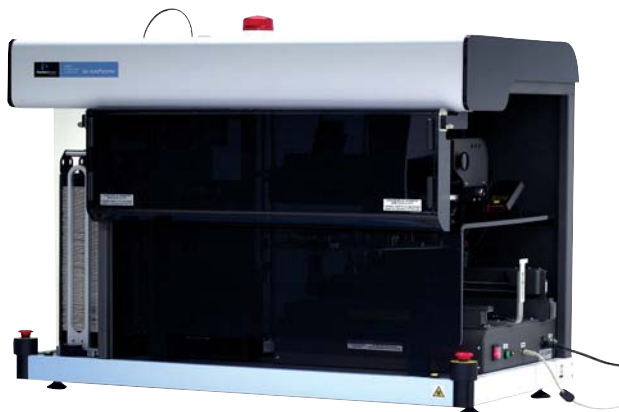


While the Wallac AutoPuncher is running no supervision is required. Nevertheless, the Workstation software clearly displays the instrument's progress.



In addition to the Workstation module, Wallac AutoPuncher also includes the independent Alert module. This is designed to attract the operator's attention whenever intervention is needed at the one or more AutoPuncher units at work in the laboratory.

The first fully automated dried sample punch for the newborn screening laboratory



1296-091 Wallac AutoPuncher

AutoPuncher instrument:	Air pump, 3.2 mm and 6.0 mm punch heads, Chute 96 (6.0), 6 Cassette Magazines, 6 Deep-well plate adapters, 6 End cards Instrument and User Manuals, Quick Guide
AutoPuncher PC:	17" flat screen, AutoPuncher software, Dell OptiPlex GX620 Stand-alone or Specimen Gate® and LifeCycle™ compatible
Calibration and Test kit	
Additional options available:	Customized bench, Cassette magazines, Deep-well plate adapters, End cards

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

©2006 PerkinElmer, Inc. All rights reserved. AutoPuncher, LifeCycle and Wallac are trademarks, and Specimen Gate 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.