



DBS PUNCHER

Punch efficiently with the DBS Puncher

The DBS Puncher is a versatile, automated instrument for punching dried blood spot samples into 96-well microtitration plates. The instrument supports a wide range

of assays employing blood spots as well as other samples dried on filter paper. The DBS Puncher is well-suited for preparing plates for feeding into an automatic immunoassay system.

Efficient punching of sample disks

The 1296-071 DBS Puncher has a user interchangeable punch head, so depending on your assay, the disks may be 1.5 mm, 3.2 mm (1/8"), 4.7 mm (3/16") or 6 mm in diameter.

Two plates may be loaded to allow simultaneous processing of different assays, and either standard or deep-well plates can be accommodated. After punching, disk detection confirms that the disk has fallen from the punch head.

The DBS Puncher is available as a standalone unit, or configured with a barcode reader and external PC with DBS Workstation software for identification of the patient sample. For process and data management, connection to an external PC also allows the transfer of worklists.

There are four selection buttons for instrument control, and the instrument is operated by hand or by foot-switch. Ease of use is enhanced by a clear double line display, and by illumination of the area being punched.

DBS Puncher

- Four different disk sizes
- Suits standard and deep well plates
- Sample identification by barcode reader
- Two plate positions for rapid handling of two assays



Suited standard or deep-well plates, and 1.5, 3.2, 4.7 or 6 mm disk sizes, the DBS Workstation option utilizes a barcode reader to identify each sample.

SPECIFICATIONS

DBS Puncher

Punch diameter: 1,5 mm (1/16")
3.2 mm (1/8")
4.7 mm (3/16")
6 mm

Plate height: 15 - 41 mm
(suits standard and deep well plates)

Operational control: 4 selection buttons for:
- operation mode selection
- plate movement
- barcode display
- operation guidance
- protocol editing

Display: 2-row, 24 character per row

Input/Output connections: RS-232 port for barcode reader
(USB with adapter)

Dimensions: Height 305 mm
Width 420 mm
Depth 480 mm
Weight approx. 20 kg

Environmental conditions: Temperature range 15-35°C
Humidity range 20-80%

ORDERING INFORMATION

1296-071	DBS Puncher
1296-8020	PC for DBS Puncher with 17" TFT Flat screen
1221-530	DBS Workstation software including barcode reader

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

Options

1296-203	3.2 mm punch head (included)
1296-2080	1.5 mm punch head
1296-204	4.7 mm punch head
1296-2070	PCR plate adapter
1296-206	6 mm punch head

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.