

## SPECIFICATIONS

### Panthera-Puncher™ 9

Newborn screening



## The new generation in dried sample punching

Panthera-Puncher™ 9 is used to punch disks from a dried sample card into the wells of up to 9 microplates. Four disk sizes are supported and two punch head sizes may be used simultaneously. Key design features include full traceability of the sample and the punching operation, and good ergonomics to support the comfort of the user.



## System Specifications

Physical dimensions	Width	876 mm (34.6")
	Depth	740 mm (29.1")
	Height	560 mm (22.0") when the instrument's feet and the monitor are set to minimum height. The monitor can be adjusted up to 100 mm higher
	Weight	45 kg (99 lb)
System	External PC, with Windows 7	
Power requirements	Power consumption	70 VA
	Mains Voltage	100-240 V, 50-60 Hz
Environmental conditions	Operating temperature	15-30 °C (59-86 °F)
	Relative humidity:	20-80 % for 3.2 mm punches and larger, 40 - 80 % for 1.5 mm punches

## Performance Specifications

### Real time visualization

- The instrument is equipped with a camera and a monitor, showing the sample card and the punching action in real-time and in color
- Easy positioning of the card
- Supports improved punching ergonomics
- The cassette/card barcode is read at the same time, with the same camera, in the same movement
- The user interface is controlled by a touch screen

### Automatically adaptable punch pattern

- Punch locations are set based on the shape and size of the blood spot
- Limit values can be set for distance from the edge of blood and distance between the punches
- After its barcode is read, the card can be punched from either side
- A fixed punch pattern may alternatively be set
- 1.5, 3.2, 4.75, 6.0 mm punch sizes are available; two punch sizes can be used in the same run

### 9 plate capacity

- 9 standard 96 well plates (ANSI SBS 1- 4), with height 14 - 16.1 mm can be loaded in the instrument at the same time. Alternatively 3 of the available plate positions may be occupied by deep well plates, with a maximum height of 44.5 mm, leaving 6 plate positions free for standard plates
- Significant improvements in the lab workflow of labs running more than 6 assays
- Punching 9 disks from one spot into different locations on 9 plates takes ~14 sec
- Minimization of static electricity by ionizer module in the punch area

### Improved ergonomics

- Designed to allow more comfortable positioning of the user's arms and shoulders
- The table may be raised or lowered electronically to suit the height of the user
- The height, distance and angle of the monitor can be adjusted.

### Simplified user maintenance

- The user can remove the punch heads from the instrument, instructions shown in the user interface
- All covers are removable to support easy access for cleaning (wiping or vacuuming)

## Safety Requirements & Conformance to Directives

The Panthera-Puncher 9 instrument meets the following safety standards:  
IEC 61010-2-101, CAN/CSA-C22.2 61010-1, UL 61010-1, EN 61326:2006, IEC 61010-1/EN 61010-1:2001.  
The punching device conforms to EU directives (2004/108/EC), (2006/95/EC) and (92/79/EEC).

## Ordering Information

### 2081-0010 Panthera-Puncher™ 9

#### including:

Punching instrument  
3.2 mm punch head  
External computer with Panthera Manager software  
External keyboard  
External mouse  
ELO touch screen (15" wide screen with stand)  
Foot pedal  
3 plate adapters  
Instrument and User manuals

#### Options:

2081-2010 1.5 mm punch head  
2081-2020 3.2 mm punch head  
2081-2030 4.75 mm punch head  
2081-2040 6.0 mm punch head  
2081-2050 Instrument table, electronic 230V  
2081-2060 Instrument table, electronic 120V

All PerkinElmer 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](http://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](http://www.perkinelmer.com/ContactUs)

Copyright ©2012, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.