

Punch intelligently



revvity

Panthera-Puncher™ 9
The new generation in dried sample punching

Panthera-Puncher™ 9



The new generation in dried sample punching

Panthera-Puncher™ 9 is the new generation dried sample punching device from Revvity. As a leader in newborn screening, Revvity has selected the technical solutions and materials needed to assure a new standard of punching excellence in terms of:

- ✓ Efficient use of sample material
- ✓ More plates simultaneously
- ✓ Enhanced traceability
- ✓ Good ergonomics

Gets your sample material to tell you more

Panthera-Puncher 9 includes a camera to provide a clear full-color view of the punching area in real-time, and an automatically adaptable punch pattern that adapts to the shape and size of the blood spot. This allows optimal use of the sample material without requiring the user to adjust the pattern to fit the area to be punched.

9 Assays from one blood card

To save time in laboratories running multiple analytes, the laboratory can prepare up to 9 microplates simultaneously. Up to 9 standard plates can be loaded at a time, and for laboratories that sometimes need to use non-standard plates, 3 of the plate positions will accept deep well plates too.

Disk sizes to suit your assay

There are 4 punch sizes, providing punched disks with diameters of 1.5, 3.2, 4.75 and 6.0 mm. Any two punch heads can be loaded simultaneously, so you can prepare all of your microplates simultaneously, even when some of the assays require a different sized disk.

Punching security through full traceability

Barcode reading takes place immediately and automatically. Samples and plates are recognized from their barcodes, contributing to loading flexibility and confirming Positive ID. The ionizer module in the punch area minimizes static electricity that could otherwise prevent the disk from falling freely, and the automatic disk recognition system then confirms that the disk has fallen into the well.

Comfortable punching

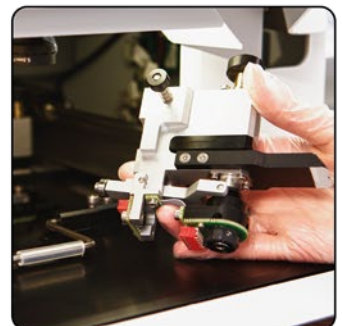
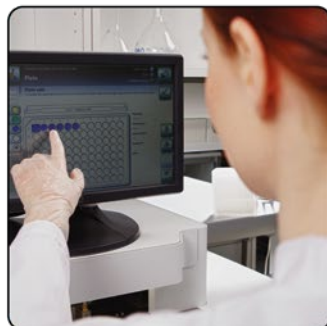
Panthera-Puncher 9 is easy to use at every stage. There are no heavy covers, and it is easy to load plates, in any order, from the front of the instrument. An electronically adjustable, ergonomic punching table is supplied as an option. Operation is by touch screen and the software has the “look and feel” of our popular Workstation software. Punching is triggered either by hand or using a foot switch.

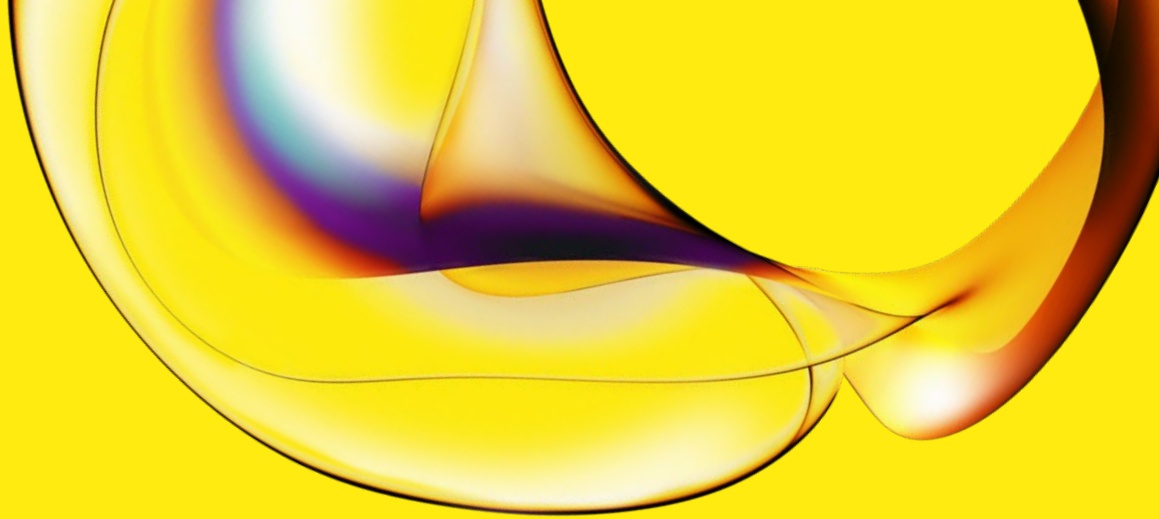
Flexible, popular design

Panthera-Puncher 9 supports a range of matrices and applications in addition to newborn dried blood spots. For some of these applications, useful features of the instrument include fixed pattern punching, punching in columns or punching blank paper.

Easy maintenance

The instrument is designed to prevent dust from accumulating and to simplify cleaning. All covers, as well as the punch head are easy to remove, with on-screen instructions and graphics to show you how.





Products may not be licensed in accordance with the laws in all countries, such as the United States and Canada. Please check with your local representative for availability.

Please note that product labeling (such as kit insert, product label, and kit box) may be different compared to the company branding. Please contact your local representative for further details.

www.revivity.com

revvity

Revvity, Inc.
940 Winter Street, Waltham, MA 02451 USA
(800) 762-4000 | www.revivity.com

For a complete listing of our global offices, visit www.revivity.com
Copyright ©2024, Revvity, Inc. All rights reserved.

1244-9975-2