

Scalable,
robust, and
automated.



revvity

EONIS™ platform



EONIS platform

Configurable SMA and SCID multiplexing from a single DBS punch

Multiplexing of SMA, SCID and XLA screening can now be done in one assay utilizing real-time PCR technology, without increasing daily hands-on work load and complexity. Designed to maximize efficiency by combining automation, multiplexing, and DNA extraction.

Following on from the launch of the worlds first commercially available (IVD) SCID newborn screening solution, Revvity continues its tradition of innovation in newborn screening with the **EONIS platform**.

The EONIS Platform has been developed by collaborating with newborn screening laboratories from around the world. The workflow is configurable to seamlessly fit in your newborn screening workflow.

Scalable, robust, and automated

- Excellent screening performance
- Ability to multiplex analytes from one punch
- Innovative analysis method calibrators
- Robust streamlined workflow with possibility for automation
- Flexibility and configurability - create the workflow that suits your needs
- Dedicated Analysis Software





Easy EONIS workflow

The EONIS assay consists of four easy steps; punching, extraction, amplification and data analysis.

Dedicated analysis software enables quantification of TREC and KREC, while SMN1 results are reported qualitatively. SMA carrier status will not be detected or reported. RPP30 is used as an internal amplification control as well as basis for the quantification. Software analysis includes automated run acceptance criteria from kit controls to ensure that the quality of measured data is not compromised.

Configurable to your throughput needs

Punching



Panthera Puncher™ 9
DBS Puncher

Extraction



EONIS DNA
Extraction kit
EONIS™ SCID-SMA kit



JANUS® Extraction
JANUS® Mini Extraction
JANUS® PCR mastermix

Amplification



QuantStudio™ Dx
Real-Time PCR

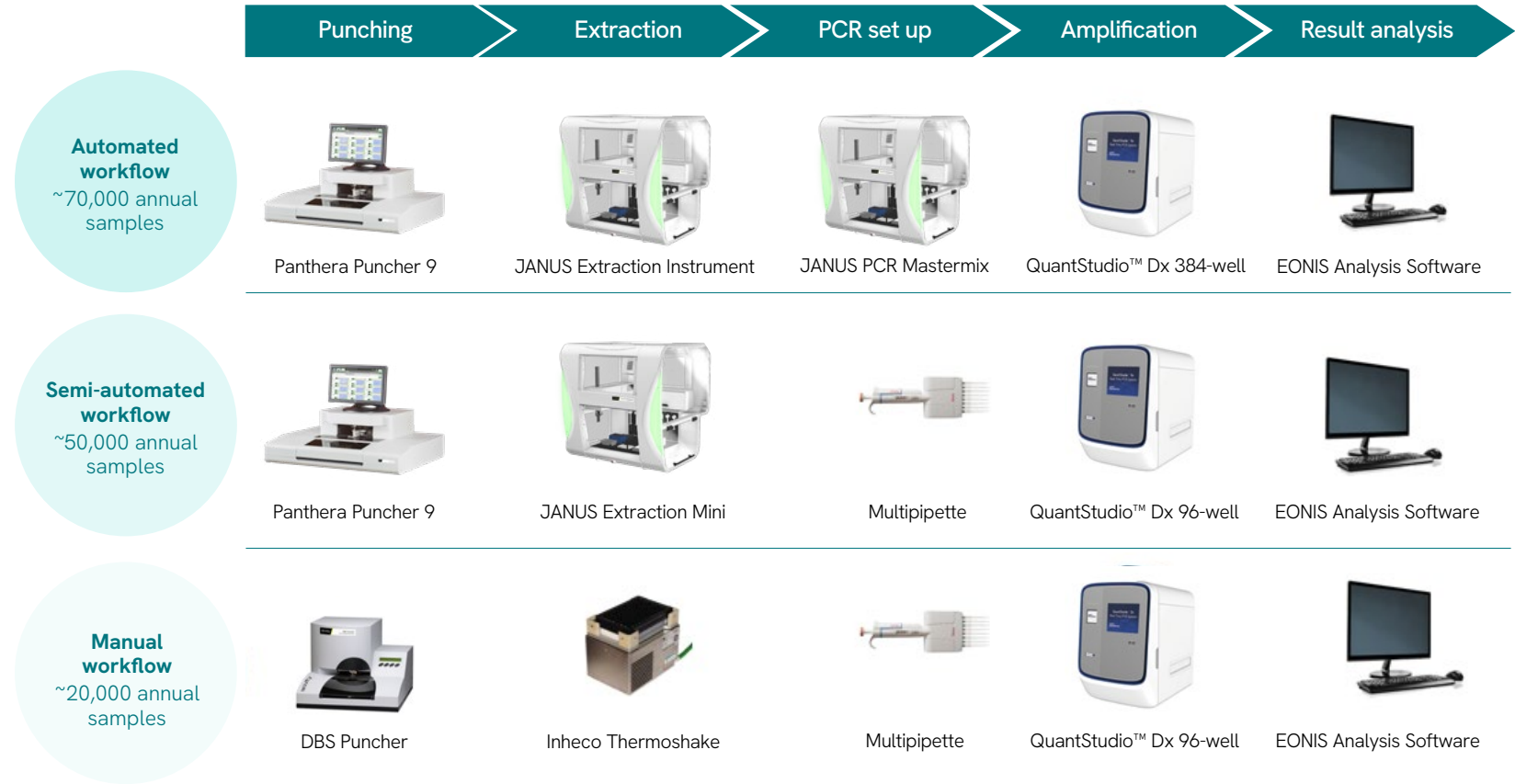
Data analysis



EONIS
analysis software



Configurable to support different throughput needs



- The design of the assay and system enables automation to further streamline
- The workflow of high-throughput laboratories without compromising the sample
- Traceability from punch to result



EONIS order guide

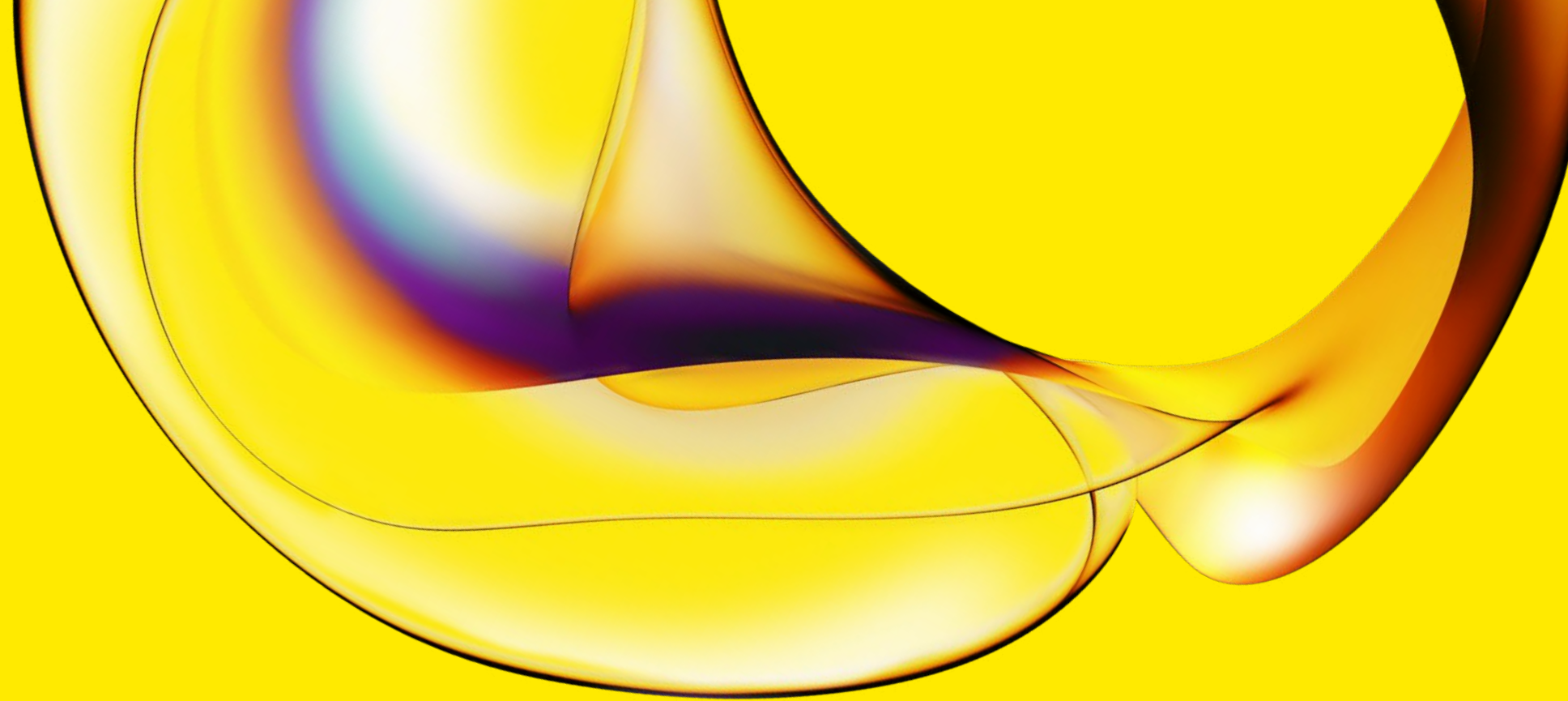
Reagents	Part number
EONIS DNA extraction kit (1152 reactions)	3240-001U
EONIS SCID-SMA kit (384 reactions)	3241-001U
EONIS SCID-SMA kit (3 x 384 reactions)	3242-001U

Software	Part number
EONIS analysis software	5020-1000
EONIS PC	5020-8010

Assay consumables	Part number
96-well PCR plate	4174-0010
384-well PCR plate	4158-0010
DNA extraction plate	4157-0010
Optical PCR seal	4159-0010
Adhesive foil seal	4156-0010

EONIS™ instruments	Part number
QuantStudio™ Dx 96 well, fast	2031-0030
QuantStudio Dx 384 well, fast	2031-0040
JANUS extraction instrument	CJM8M03
JANUS Mini extraction instrument	CJSM003
JANUS PCR mastermix instrument	CJS8003

Calibration consumables	Part number
96-well calibration plates kit	3243-0010
96-well calibration plate, background	3248-0010
96-well spectral calibration plate with Cy5 dye	4172-0010
96-well spectral calibration plate with Cy5.5 dye	4173-0010
384-well calibration plates Kit	3225-0010
384-well calibration plate, Background	3226-0010
384-well spectral calibration plate with Cy5 dye	4178-0010
384-well spectral calibration plate with Cy5.5 dye	4179-0010



www.revivity.com

revivity

Revvity, Inc.
940 Winter Street, Waltham, MA 02451 USA
www.revivity.com

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

1254004

For *in vitro* diagnostic use.
Products may not be licensed in accordance with the laws in all countries.
Please check with your local representative for availability.