

Faster molecular
analysis made
easy, for all.



revvity

EONIS™ Q qPCR Workflow
Screen for SCID and SMA

Simplicity and speed, streamlined

Initiatives to screen for SCID (severe combined immunodeficiency) and SMA (spinal muscular atrophy) are gathering huge pace across the globe, yet challenges exist in overlaying new solutions into very busy, crowded newborn screening labs.

For both inherited conditions, immediate detection is critical to advancing a positive outcome.

For SMA, disease modifying therapies exist to stop progression of disease,¹ and for SCID,

immunoglobulin treatments combined with stem cell therapies can potentially cure a child, if intervention arrives in time.²

The EONIS Q system simplifies and streamlines molecular testing for SMA and SCID with an innovative workflow, comprising the EONIS™ Q96 instrument, the EONIS™ SCID-SMA kit and dedicated EONIS™ EASI™ software.

Three steps for working smarter, not harder

With just three steps, workflow complexity is reduced, minimizing the need for specialist skills sets, while gaining fast and accurate results of controls and samples.

Turnaround time for the EONIS Q workflow, from sample-in to result, is approximately three hours. In comparison to conventional wet qPCR solutions this can offer valuable time saving benefits.

1. Punch samples

Single punch works for up to three analytes.



2. Elute

No wash steps required, saving you valuable time.



3. Analyze

Fast results of controls and samples.



How we've simplified the workflow

- ✓ Reagents are dried on the plate: no need for a master mix preparation
- ✓ Multiplex from a single dried blood spot punch
- ✓ Quantification without calibrators
- ✓ Analyze over 500 samples per day

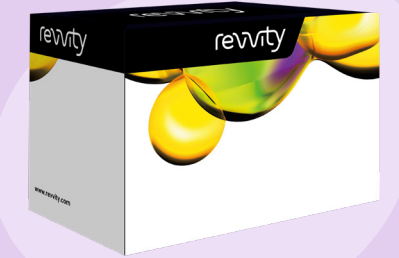
The EONIS Q system

System simplicity to maximize your runs!

The EONIS™ Q is designed to be compact, offering a streamlined workflow with a swift 3-hour turnaround, eliminating wash steps and reducing the need for extensive pipetting and centrifuging processes.

EONIS Q reagents are pre-plated and dried to limit any chance of contamination, meaning no pre-PCR clean room needed.

EONIS Q kits include lot specific calibration information included in the pack insert allowing for more tests to be run at one time.



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.

www.revity.com

EASI Software

Uniquely customised for newborn screening laboratories

- Easy to use without any previous experience of qPCR software
- The same software is used for managing the PCR run and analysing the results

Seamless integration

- Connects to LIMS system easily with no extra tools needed to run analysis

Time to report

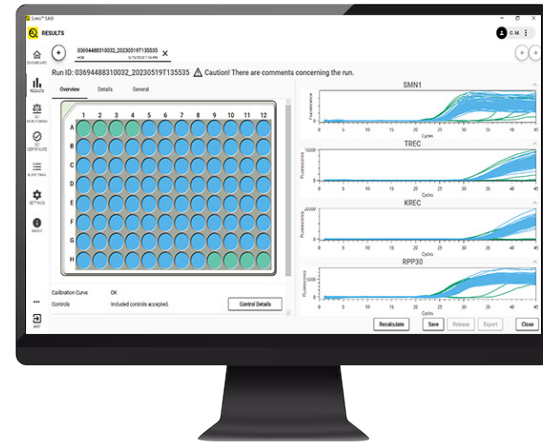
- Fast reporting: no manual data handling that requires extra time

Troubleshooting packages

- With just one click, you can easily share files with our customer support if you need assistance

Quality control monitoring

- Built in QC monitoring ensures you can track your QC levels over time: no need to keep any separate files for tracking



EONIS Q screens for diseases that can potentially be treated or cured.



Spinal muscular atrophy (SMA)	Severe combined immuno-deficiency (SCID)	X-linked Agamma-Globulinemia (XLA)
SMN1	TREC	KREC
1 : 7,000 - 10,000	1 : 50,000	1 : 100,000 - 200,000
Modulating SMN2 Gene therapy	Hematopoietic stem cell transplantation, bone marrow transplantation, gene therapy (ADA-SCID)	Intravenous immunoglobulin therapy, prophylactic antibiotics



The EONIS Q system and workflow has been designed to make screening for SCID and SMA easier by optimizing time, space, cost and training!



Enhances the whole lifetime cost of the system

Uses standard pipettes, it all fits on a small table



Decreases space needed

No need for a clean pre-PCR space



Minimizes hands-on time

Compared to wet chemistry workflow



A simple system improves potential training times

Making staff training easier



Improve efficiency with unconfigured workflow checks



Removes the need for constant data transferring checks

The system seamlessly connects with the LIMS system

Screening for SCID and SMA has now been made easier for all labs

Start your SCID and SMA screening program today by contacting your local Revvity representative.

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.

Ordering information

Instruments

Product name	Product number
Eonis Q96 instrument	2044-0010
Eonis EASI software	2044-3010
Eonis Q96 Dx EASI software licence	2044-3020
Eonis Q PC, monitor and barcode reader	2044-8010
96-well optical test plate (for installation and PM only)	X04-10-310P-105-14
TriNEST microplate incubator shaker	1296-0050



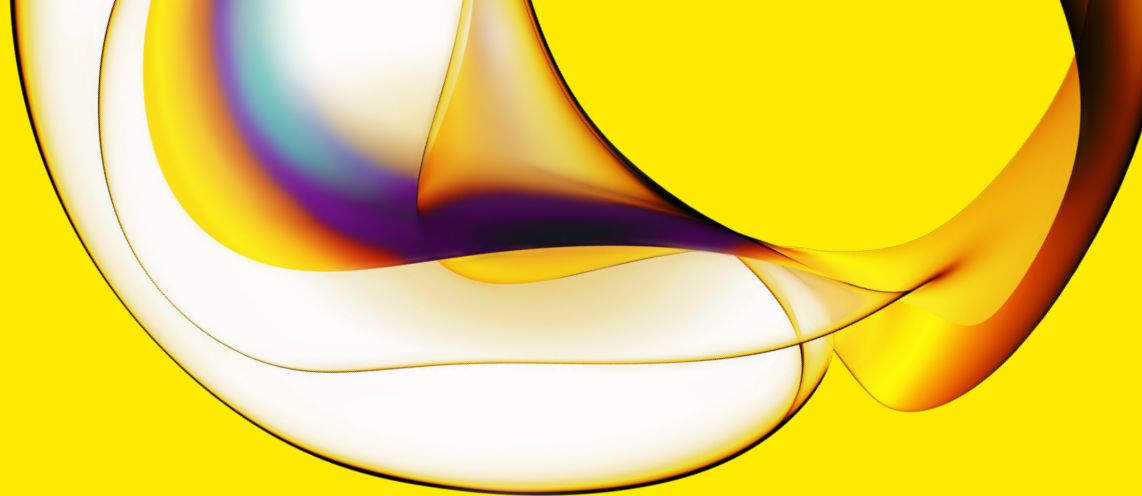
Reagents

Product name	Description	Product number
Eonis SCID-SMA kit	4x PCR plates; 96 reactions/plate, 88 samples/plate 1x bottle Elution Solution 2x Kit controls 1x Zero control	3241-0020

Consumables

Product name	Description	Product number
DNA elution plates	20 plates	4183-0010
Adhesive foil seal	100 plates	4156-0010
Optical PCR seal	100 plates	4197-0010

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.



- 1 *Spinal muscular atrophy: Screen at birth, save lives.*
Whitepaper Version 2, 25 November 2021 European Alliance
for Newborn Screening in SMA
- 2 *Classic Severe Combined Immunodeficiency* Rebecca Buckley,
MD, Duke University School of Medicine, Durham, North
Carolina, USA Jennifer Heimall, MD, Children's Hospital
of Philadelphia, Philadelphia, Pennsylvania, USA, Immune
Deficiency Foundation Patient & Family Handbook For Primary
Immunodeficiency Diseases 6th Edition

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

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 ©2023, Revvity, Inc. All rights reserved.

1151724