

Step into newborn sequencing.

Powering the next wave of genomic research.



Newborn research stands at the dawn of a whole new era. To date, standard research efforts have documented hundreds of disorders in the context of newborn health. However, less than 80 of them are actively investigated. Now, with the advent of sequencing technology, the limitations have been lifted.

Bringing simplicity to sequencing and featuring a panel of 390+ genes, Revvity's newborn sequencing workflow is powering the next frontier of research.

Unleash your laboratory's potential. Explore Revvity's newborn sequencing research workflow today.

The human genome at a glance:

3 billion
DNA base pairs

~20,000
genes total*

6 feet tall
human DNA
stretched out flat



*Estimated number of genes that produce proteins

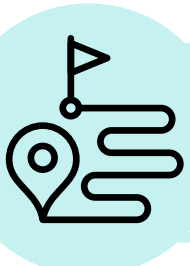
Newborn sequencing with Revvity

A solution for every step



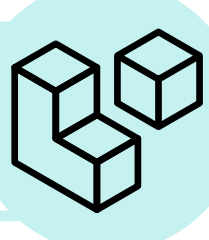
From dried blood spot card to final data analysis.

Revvity's lab workflow



End-to-end solution

Customizable workflow



Hands-on NGS expertise

Multiple starting points



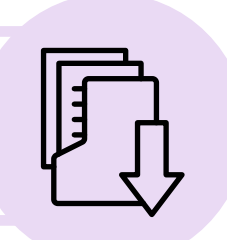
Adopt full solution (DIY)

End-to-end service



Research report service

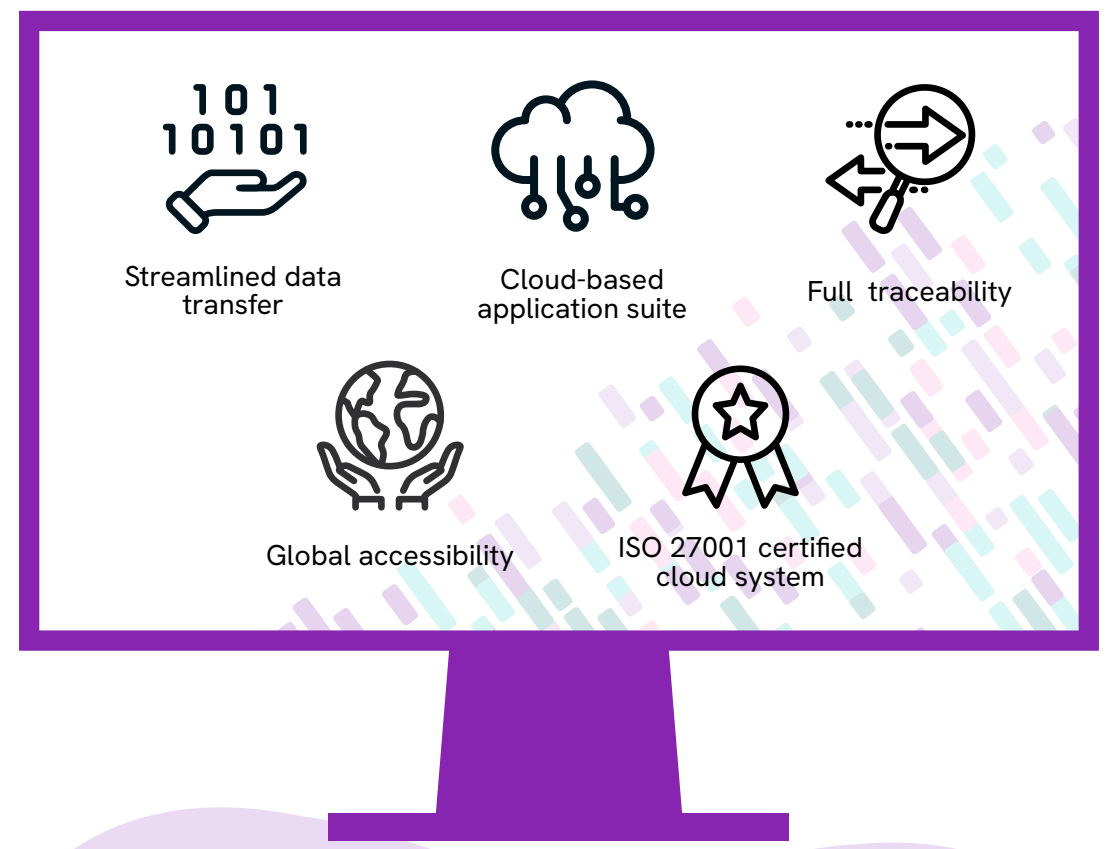
Data generation service



Complete software suite

- LIMS
- Analysis
- Interpret

Software designed for the lab



Key specifications

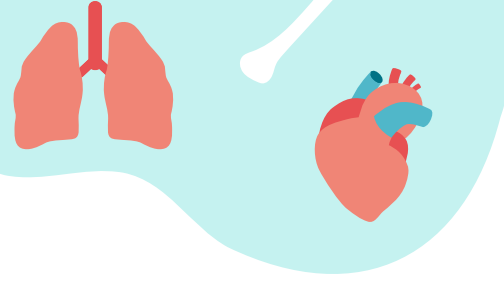
- Software engineered by laboratory specialists
- Compatible with diverse laboratory setups
- Runs on dried blood spot samples, whole blood, buccal swaps and saliva

Analysis and interpretation

>390:
number of genes we test for



17
research areas



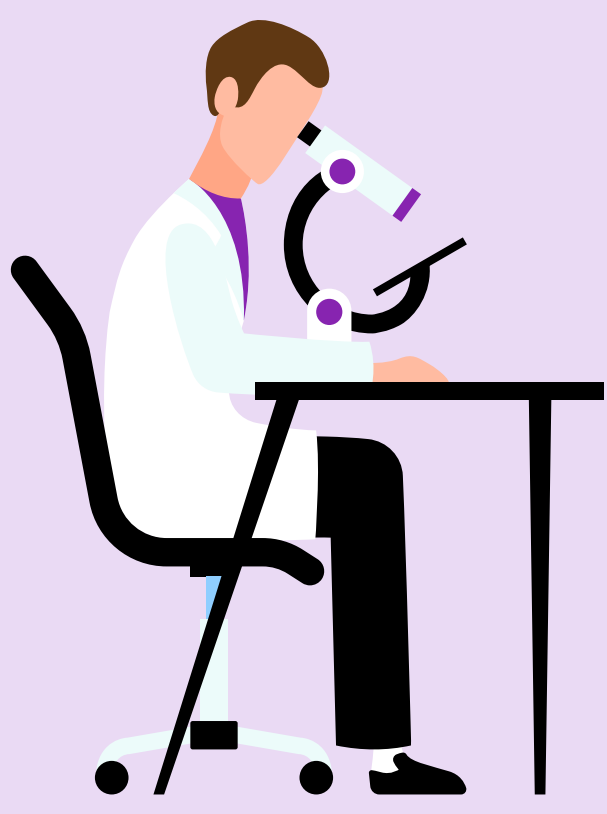
Carefully curated panel



>87K :
pre-curated variants
with written interpretation



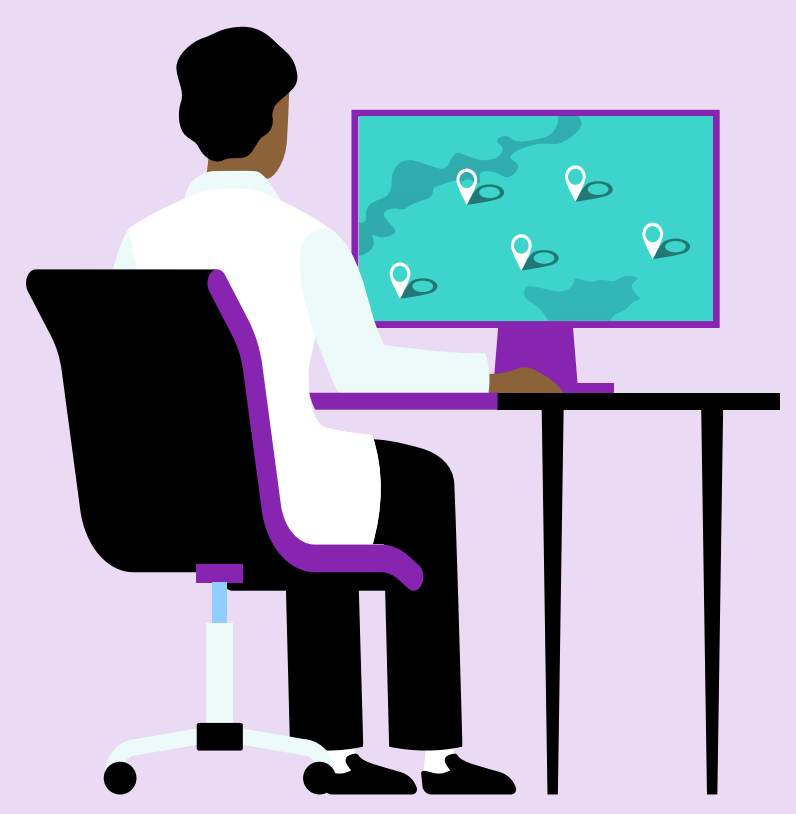
Revvity lab service



Over **40K** samples processed



Validated by **CLIA, CAP, NY State, ISO15189** standards



5 Revvity NGS labs globally

Why Revvity?



> 40 years experience, the leaders in newborn screening



> 75 years providing dynamic research lab solutions



190: number of countries empowering academic & translational research

See how your lab can step into sequencing.

[Learn more](#)

For research use only. Not for use in diagnostic procedures.

revvity

Copyright ©2024, Revvity. All rights reserved. Revvity is a registered trademark of Revvity.

1466204