

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, Revvity's newborn sequencing workflow features a robust gene panel curated by genomic experts, powering the next frontier of research.

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

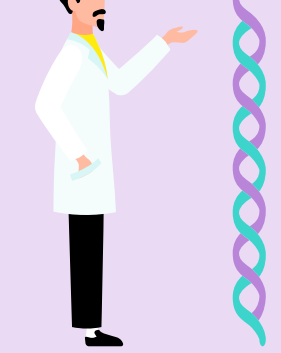
[Learn more](#)

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

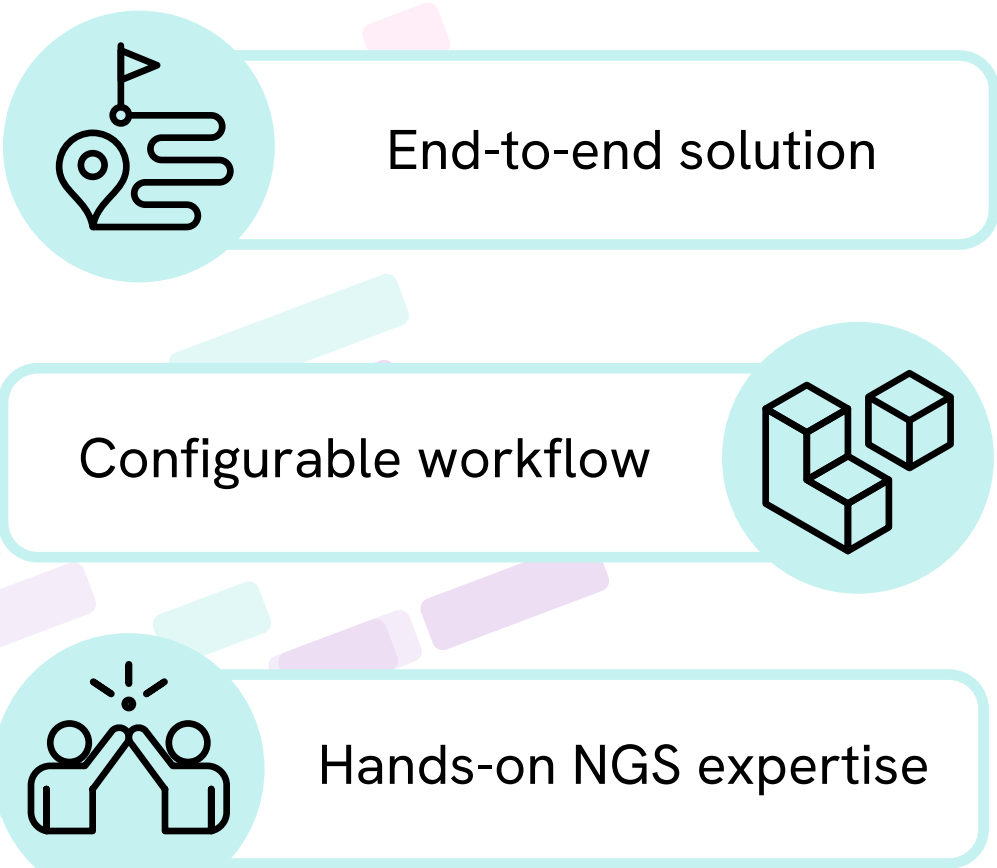
Newborn sequencing with Revvity

A solution for every step



From dried blood spot card to final data analysis.

Revvity's lab workflow



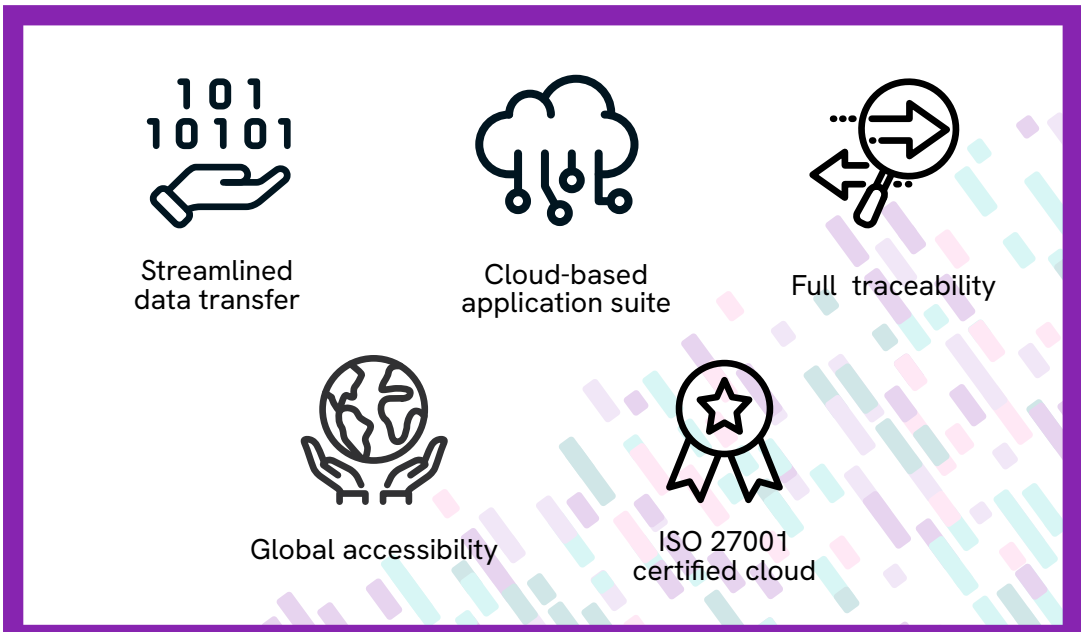
Multiple starting points



Complete software suite

- LIMS
- Analysis
- Interpret

Software designed for the lab



Key specifications

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

Analysis and interpretation



Revvity lab service



Over **40K** samples processed

Validated in accordance with **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.

*This testing service has not been cleared or approved by the U.S. Food and Drug Administration. Testing services may not be licensed in accordance with the laws in all countries. Licenses, certifications, and specific test offerings vary by laboratory location. This content is provided for informational purposes only, not as medical advice. It is not intended to substitute the consultation, diagnosis, and/or treatment provided by a qualified licensed physician or other medical professional.

revvity