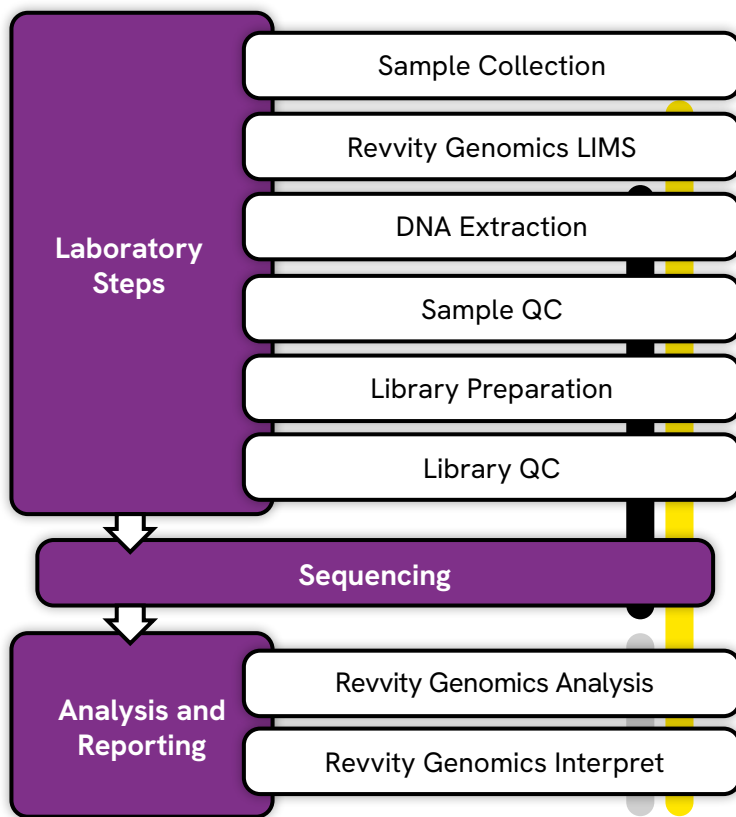


Discover the future of newborn sequencing research

A customizable, end-to-end sequencing workflow equipped to your laboratory's needs.



Newborn research, reimagined.

Across the globe, there are technological gaps which can limit the scope of newborn sequencing research. With decades of laboratory expertise, Revvity's end to end sequencing workflow - from blood spot card through final analysis - bridges these gaps, empowering newborn research beyond expectation.

Note: Workflow is compatible with a host of lab instruments to best suit your needs. Please reach out to your Revvity representative for more information.

Options for Outsourcing

- Full end-to-end service
- NGS data generation service or flowcell rental
- Reporting service

Power your newborn sequencing research with Revvity today.

Learn More

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

revvity