



Discover the future of newborn screening research.

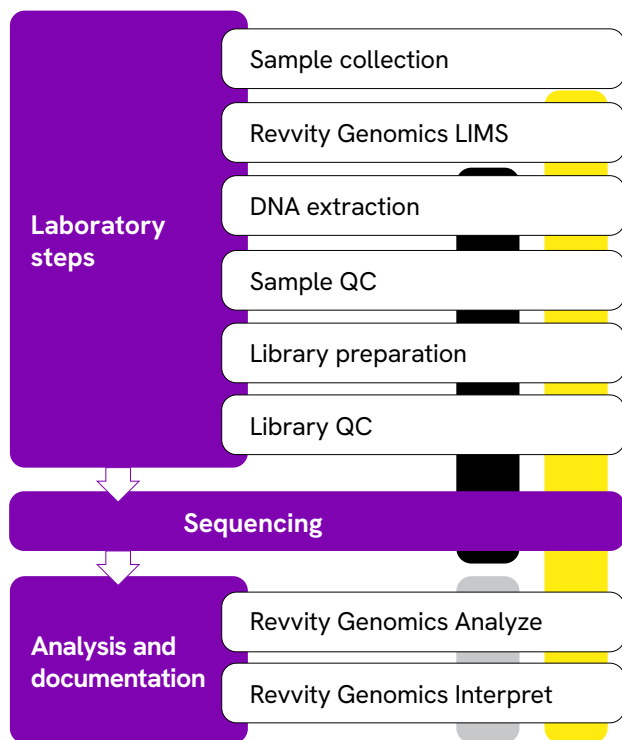
One configurable, end-to-end sequencing workflow equipped to your laboratory's needs.

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.



Options for outsourcing:

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



Scan or click
for more information



Scan or click
for more information

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.