



Is your laboratory ready for newborn sequencing?

We're here to help. Use our NGS readiness checklist to evaluate your sequencing strategy, operational capacity, and infrastructure. Identify strengths, gaps, and the most efficient route to implementing or expanding newborn genomic research.

Criteria	Consideration	Laboratory notes
Foundational elements	Project objectives: pilot study etc.	
	Genomic scope: WGS, WES, targeted panel	
	Annual sample volume	
Infrastructure	Sample collection and transport	
	Integration with existing LIMS and/or instrumentation	
	Lab environment: temperature-controlled storage, clean room environments, physical space requirements	
Instrumentation	Sequencing platforms	
	Sample preparation automation (puncher, DNA extraction, library prep)	
	Supporting equipment (QC instruments)	
Software	Genomics LIMS	
	Bioinformatics data analysis	
	Variant interpretation, classification and reporting	
	Infrastructure and data management	
Expertise	Regulatory compliance, security and privacy	
	Laboratory tech training	
	Clinical genetics knowledge	
	Bioinformatics expertise	
Investment	Data management specialists	
	Capital equipment	
	Genomic software	
Format	Data storage costs	
	In-house workflow installment	
	Send-out lab services	
	Development of a lab-in-a-lab	

Questions about your lab's path to NGS? Connect with our newborn genomic experts to discuss sequencing strategies, infrastructure planning, informatics and more.

Contact us

