



**Agilent**

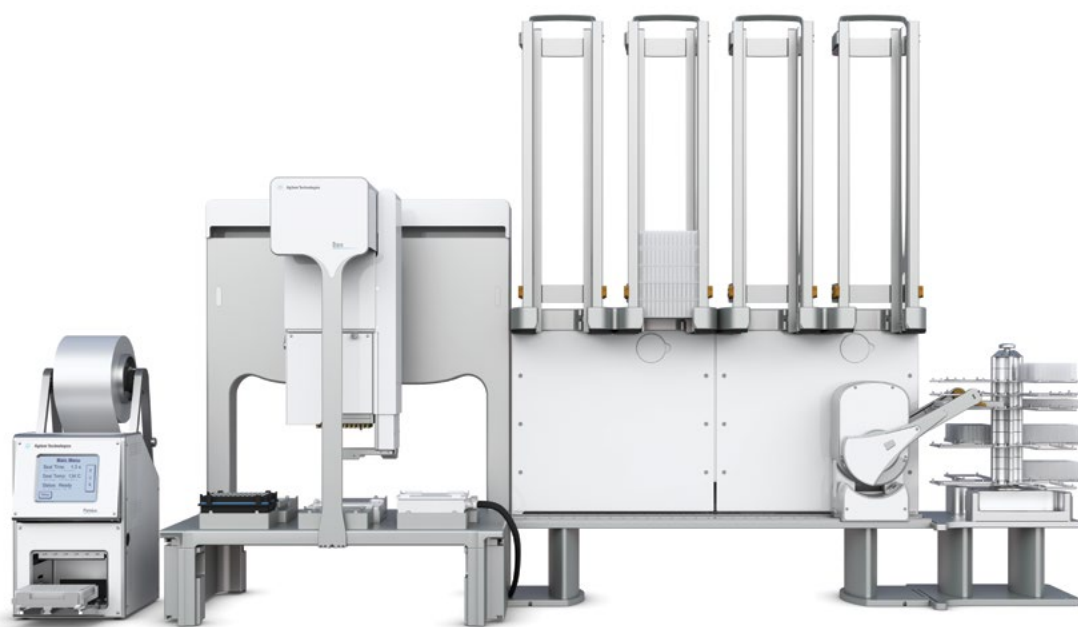
Trusted Answers



# Accelerate Your Research

## Agilent Automated NGS Sample Prep

- **Increase walk-away time:** Benefit from ready-to-use automation protocols.
- **Maintain consistent data quality:** Perform pipetting steps with high precision, every time.
- **Adapt to changing needs:** Effortlessly run multiple workflows on the same platform and scale up throughput with ease.



For Research Use Only. Not for use in diagnostic procedures