

Quanterix™

The Science of Precision Health

Quality Policy

Quanterix is dedicated to the sustainable delivery of high quality, innovative biomarker measurement products and testing services that meet our customer needs and regulatory requirements.

Sustainable quality is accomplished via maintenance of an effective Quality Management System that includes the following:

- Executive commitment to an environment of operational excellence, continuous improvement and customer satisfaction
- Recruitment and retention of the best and brightest workforce that embodies the principles of Accountability, Teamwork, and Trust
- Rigorous management of our product design and development process
- Establishment of appropriate goals throughout the organization to achieve the quality objectives
- Continuous monitoring and improvement of our process metrics, product performance and quality objectives
- Partnerships with our suppliers and customers for feedback to achieve product and service excellence
- Supporting engagement of our workforce throughout the Quality Management process

*QUANTERIX Sr. Management Team has developed this quality policy. The policy is explained and discussed at the general orientation training given to all existing and new employees. The policy is also posted in locations throughout the company.

Members of the Sr. Management Team are part of the Management Review process, including reviews of the Quality Policy for adequacy, effectiveness, and relevance