

## QIAGEN Digital Insights



Unlocking the secrets of 'omics data for better outcomes.

---

**'Omics, the vast amount of raw data generated by current molecular diagnostics, such as next-generation-sequencing, is colossal, complex and changing – how can this change translate into results ? QIAGEN Digital Insights (QDI) is the leading provider of genomic and clinical knowledge, analysis and interpretation tools and services for scientists and clinicians. QDI integrates the QIAGEN acquisitions of Ingenuity, CLC, OmicSoft, BIOBASE and most recently N-of-One under the QIAGEN brand. QDI offers the largest publicly available source of expert curated molecular information in the world to accelerate Sample to Insight.**

---

Bioinformatics data is complex and expanding at an unprecedented rate. Estimates indicate there are over 28,000 peer reviewed journals with 2.5 million new scientific publications a year growing at over 8-9% per year. Staying up-to-date and informed is a challenge for every scientist and clinician. Digital Insights uses over 350+ experts and advanced AI software to develop the QIAGEN Knowledge Base. Supporting exploratory, analysis and reporting software and services, the Knowledge Base is foundational in enabling scientists and clinicians to see new connections, improve understanding and advance interpretations.

### QIAGEN sets standards in bioinformatics

Digital Insights is applying its products, knowledge and services to provide Sample to Insight offerings as well as platform agnostic solutions to help transform the workflows and productivity of researchers and clinicians. Discovery products support a wide range of research workflows from gene expression analysis to bacteria and virus characterization. While combining QIAGEN's new QIAseq Multimodal Panels for simultaneously sequencing of RNA and DNA from a single sample, QIAGEN Digital Insights software can rapidly analyze raw NGS data, filter and identify variants

before delivering expert curated variant interpretation reports leveraging the Knowledge Base. The result is customers have seen Sample to Insight turnaround times up to 85% faster than other methods.

QIAGEN Digital Insights brings together over 90,000 product users and 3 product portfolios to meet the growing demand for molecular insights; QIAGEN Discovery Insights, QIAGEN Clinical Insights and QIAGEN Real-World Insights. Together these portfolios support bench to precision medicine for insights from early discovery, to clinical trial optimization and genetic variant interpretation for hereditary and oncology. To date QIAGEN Digital Insights clinical products have been used in over one million pathway studies and 1.7 Million hereditary and oncology molecular profiling cases to make improvements in life possible.

### QDI proven depth and breadth of Insights

Data is one of the few commodities that the more people you share it with people the more valuable it becomes. Digital Insights is focused on building our Knowledge Base to deliver a single source of trusted knowledge curated from millions of scientific findings and patient studies that is easily accessible to everyone. QDI customers can access information to improve insights and accelerate decisions from over 40 expertly curated databases, covering 20 million findings and 33,000 diseases to improve scientific and clinical outcomes. It's why today we have over [37,000 references](#) that publicly cite Digital Insights products and services.

Examples of QIAGEN's Digital Insights areas of application:

- **QIAGEN Discovery Insights:** Research and discovery – build on the knowledge of millions to unlock new scientific discoveries
- **QIAGEN Real-World Insights:** Translational research and clinical development – identify and prioritize target populations for precision therapeutics
- **QIAGEN Clinical Insights:** Public health, testing laboratories – empower clinical genomics labs with trusted insights for precision medicine

