

QIAGEN 2016: Results and Important Events

	2016	2015
Net Sales (US\$M)	1,34 (+6% CER*)	1,281,2
Adjusted EPS** (US\$)	\$1.14	\$1.05

Dec 12	QIAGEN welcomes new evidence-based tuberculosis testing guidelines from U.S. professional societies: New guidelines recommend broader use of modern blood TB tests like QuantiFERON-TB Gold over the 100-year-old tuberculin skin test
Dec 1	QIAGEN launches unique CALR mutation assay for MPN diagnosis
Nov 9	QIAGEN GeneReader NGS System to relaunch in U.S. market
Oct 20	QIAGEN launches QIAcscout for single-cell isolation
Sep 29	QIAGEN and ICMP to develop first NGS workflow for identification of missing persons: Collaboration targets "Sample to Insight" sequencing solutions to enhance research and applications in human identification
Sep 8	QIAGEN welcomes U.S. guidance for latent TB testing in primary care
July 27	QIAGEN launches novel integrated system for collection, stabilization, transport, storage and isolation of circulating cell-free DNA
July 27	QIAGEN launches "Digital NGS" solutions with unique Molecular Index Technology to reliably detect all genetic variation
June 2	QIAGEN forms Lung Cancer Expert Alliance: Key opinion leaders' expertise to guide design of new gene panels for GeneReader NGS
June 1	QIAGEN launches liquid biopsy workflow for GeneReader NGS System: Actionable Insights Tumor Panel enables monitoring of cancer biomarkers from plasma samples
May 25	QIAGEN partners to develop predictive tests in oncology: Collaboration with Therawis addresses unmet need to guide clinical decisions in breast cancer
April 13	QIAGEN partners to develop novel companion diagnostic in lung cancer: Collaboration with Mirati Therapeutics targets MET gene mutations to personalize treatment
April 4	QIAGEN's GeneReader NGS System wins international design distinction
Jan 11	QIAGEN reaches milestone with portfolio of master collaboration agreements for pharma companion diagnostics

* CER, Constant Exchange Rates **For further information on adjustments, please refer to the reconciliation tables in QIAGEN's Q4 2016