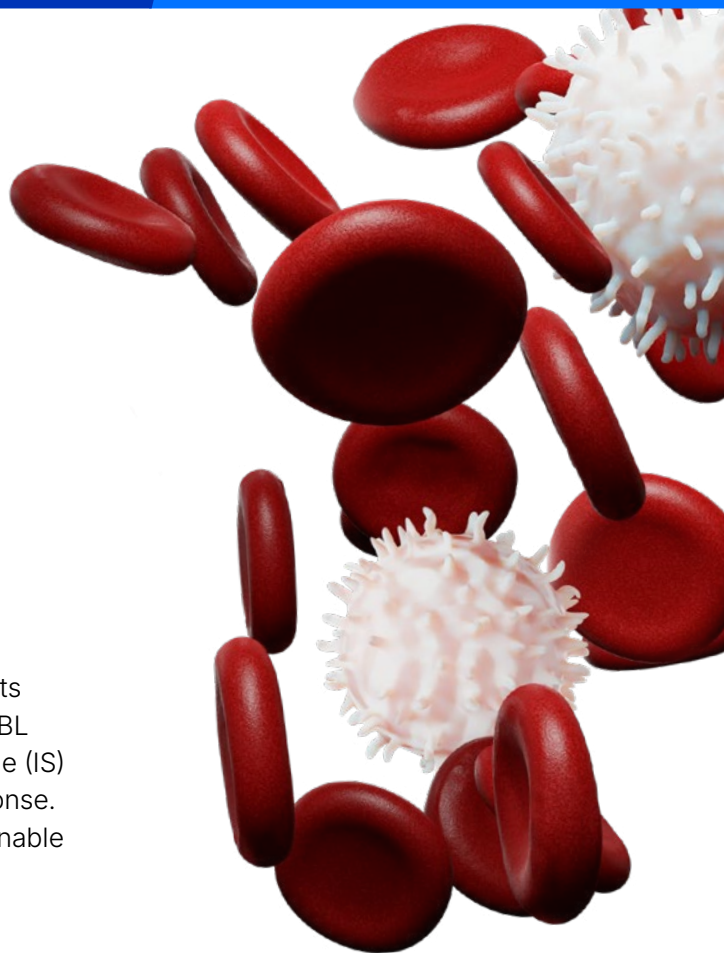


biotechne
diagnostics™

BCR-ABL Monitoring in CML Patients *with Results Aligned to the International Scale*

The QuantideX® qPCR BCR-ABL IS Kit provides confident results as an IVDR Class C-certified CML monitoring assay for BCR-ABL major breakpoints (e13a2, e14a2). With direct International Scale (IS) reporting, laboratories can easily assess deep molecular response. The multiplex assay design and integrated analysis software enable fast, accurate reporting in about four hours.



Reduced Complexity

- Direct reporting on the IS with four calibrators traceable to the WHO primary standards
- All-in-one kit and software from a single vendor

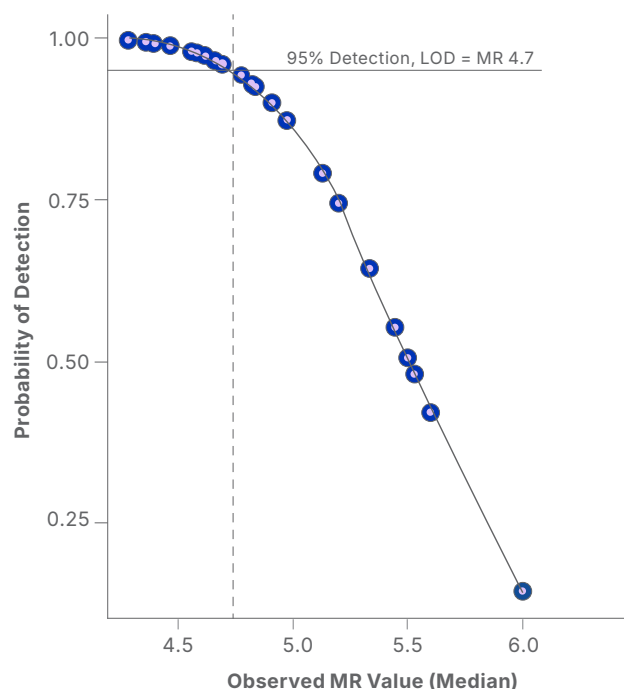
Optimized Workflow

- Multiplex design detects BCR-ABL and ABL in the same reaction, with ability to run up to 49 samples per plate
- Simplified reporting with included QuantideX Reporter software

Quality Results

- Limit of Detection of MR4.7 (0.002% IS) was confirmed in clinical human blood samples, not cell lines
- Trusted, reproducible results with Class C IVDR certification

FIGURE 1.



MR4.7 with 95% detection was determined from 1,678 valid data points from clinical samples, from 40 test runs by four operators over two kit lots, 15 days, and four instruments.*

*Brown JT, et al. J Mol Diagn. 2019 Jul;21(4):718-733.

Streamlined and Scalable Workflow

FIGURE 2.

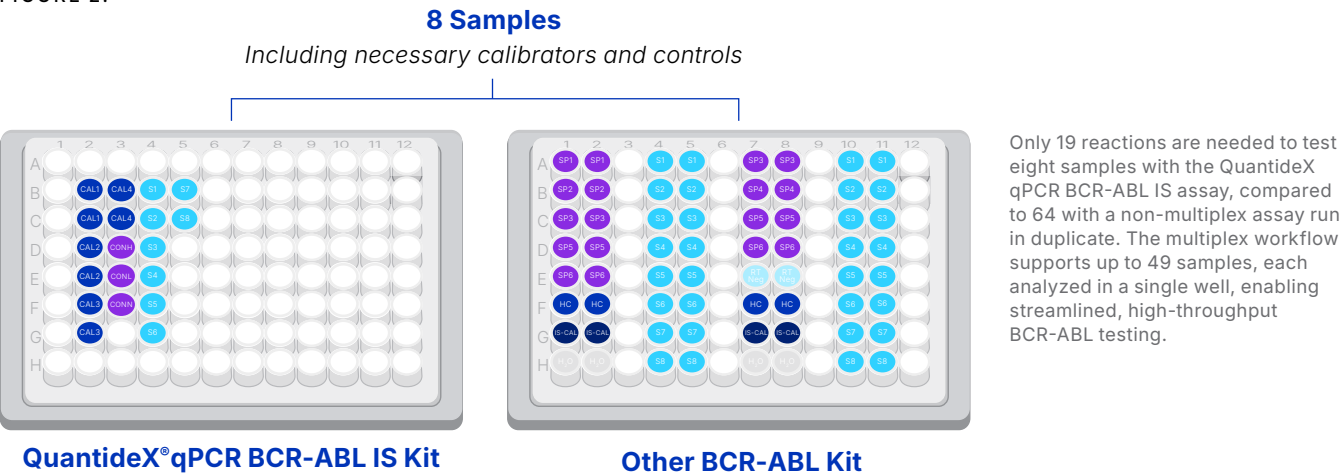
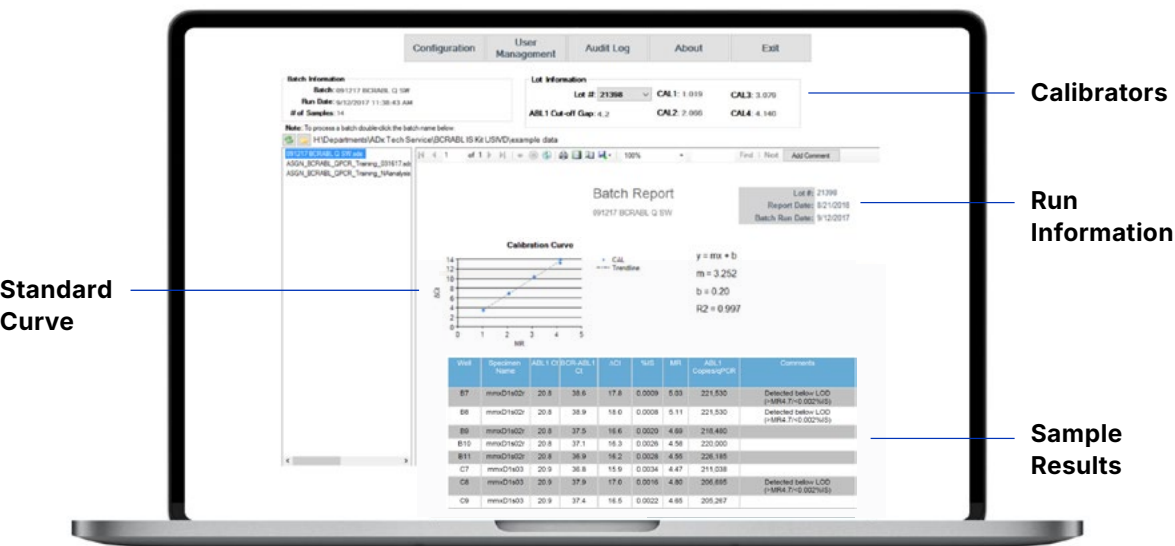


FIGURE 3.



The QuantideX Reporter automates qPCR BCR-ABL IS result reporting, performing all calculations and generating batch reports with the Standard Curve, MR values, and %IS for controls and patient samples. Reports, including the Standard Curve, are produced in PDF, XML, and CSV formats for easy LIMS integration.

Ordering Information

Part Number	Product Description	Number of Reactions
86003	QuantideX qPCR BCR-ABL IS Kit (IVDR-certified)	60



Contact Us
[bio-technediagnosics.com](https://www.bio-technediagnosics.com)
Scan the QR code

biotechne diagnostics

CE-IVD for US Export Only.
Bio-Techne Diagnostics™ and Bio-Techne® are trademarks or registered trademarks of Bio-Techne Corporation and affiliated entities. All other trademarks, service marks, and trade names are the property of their respective owners. Any use of third-party names, logos, or marks does not imply affiliation, sponsorship, or endorsement. © 2026 Bio-Techne.