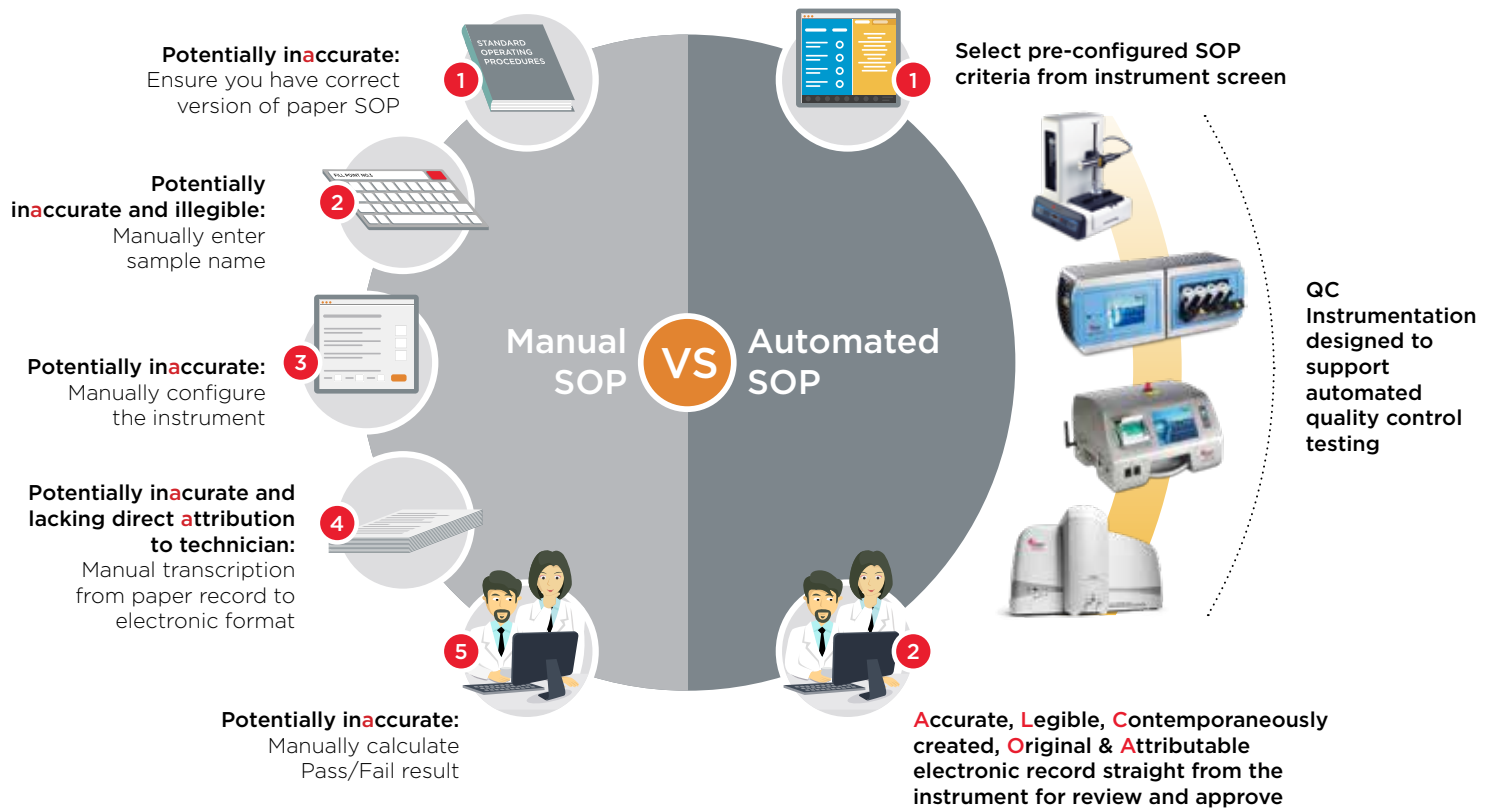


Automating Quality Control Testing for Improved 21 CFR Part 11 and ALCOA Data Integrity



How Beckman Coulter Life Sciences QC Suite instruments help you comply with 21CFR part 11 ALCOA

- **Attributable** - we offer electronic signatures, so that the record is attributable to the technician
- **Legible** - electronic record straight from the instrument avoids illegible manual writing and typos
- **Contemporaneously** - avoid the common practice of collating multiple paper results over a period of time then typing them into an electronic record at a later date, giving rise to opportunities to lose pieces of paper, or forgetting which result relates to which measurement. Our electronic record straight from the instrument allows you to create the electronic record in the same timeframe that the test is done, i.e. contemporaneously
- **Original** - avoid the common industry practice of starting with paper print-outs and then manually typing them in. Our instruments allow you to create the electronic 'original' straight from the instrument
- **Accurate** - by avoiding manual data entry and automatically generating a Pass/Fail report according to your SOP criteria in electronic format our instruments help you to ensure accuracy of the final record

Discover more at beckman.com/alcoa



© 2019 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at beckman.com

PART-5453IFG06.19