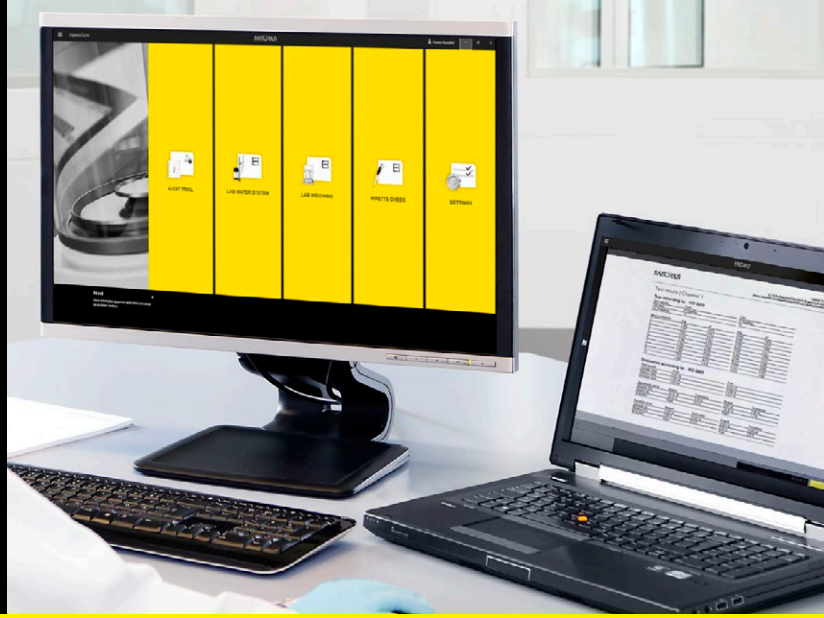


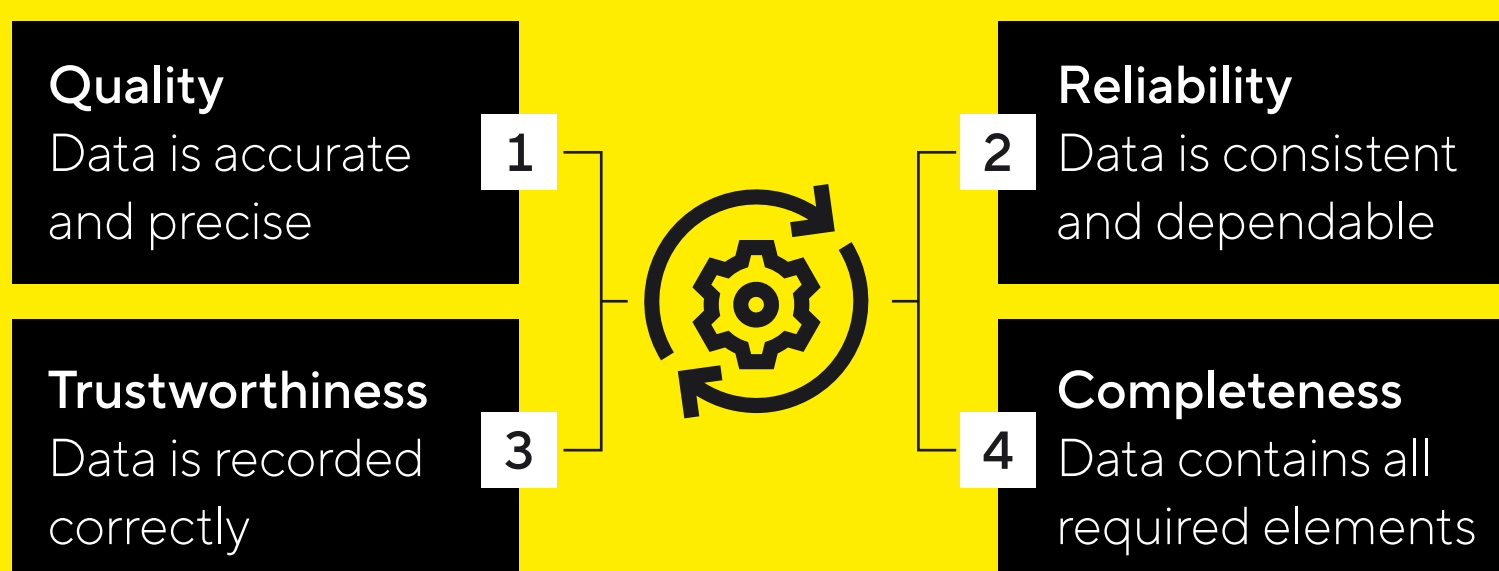
# The Path to Rock-Solid Data Integrity

Building Credibility into Scientific Research and Pharmaceutical Manufacturing

Data integrity is crucial for achieving reliable outcomes. It supports regulatory compliance, product quality, and ultimately, patient safety.



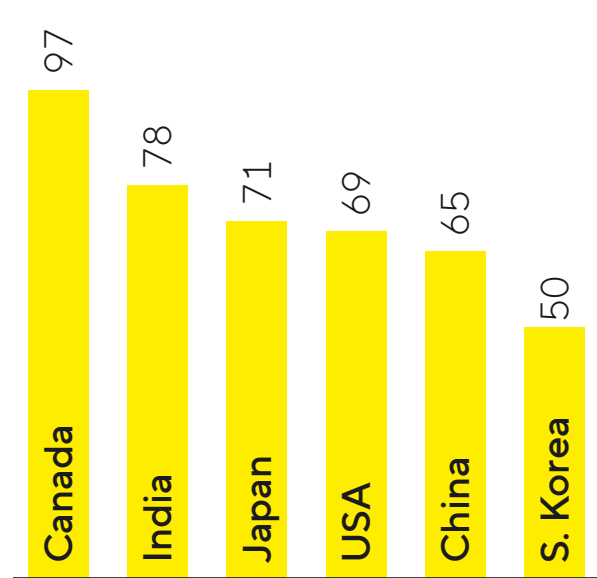
## What is Data Integrity?



## FDA Targets Data Integrity Violations

- Data loss or manipulation
- Improper batch documentation
- Insufficient audit trail
- Procedure noncompliance
- Incomplete data transfer
- Loose paper data

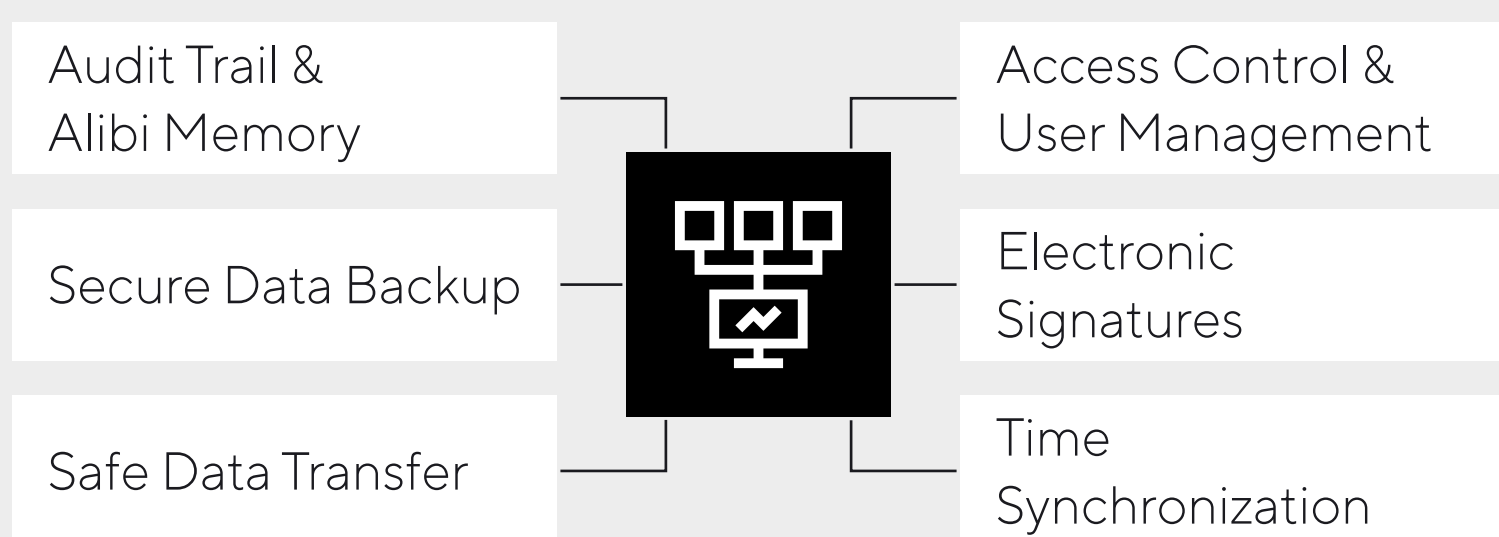
\* [Chart data published in 2021 article](#)



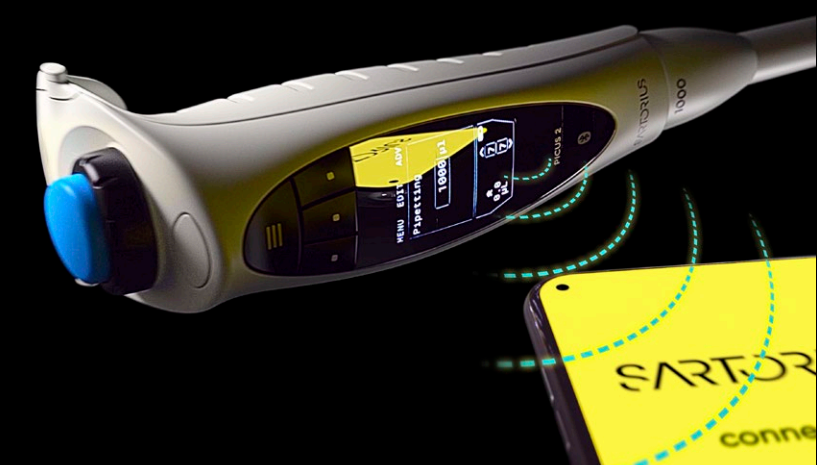
% Drug GMP Warning Letters Citing Data Integrity Violations

## Automating Compliance

Modern lab devices simplify data integrity compliance by replacing manual, paper-based procedures with automated technical controls.



## Connected Solutions



### Electronic Picus® 2 Pipettes

- Open connectivity
- Bluetooth-enabled
- Password protection
- Calibration reminders
- Pipetting app



### Cubis® II Lab Balances

- 21 CFR Part 11, EU Annex 11 compliance
- Complete traceability
- Submission-ready data
- User management controls
- SOP integration

## Ingenix Suite Digital Fleet Management



- External balance calibrations and weight set management
- Open software solution
- ALCOA+ principles; 21 CFR 11 and EMA Annex 11 compliance

- Pipette performance checks and calibrations

**Find out more:**  
Ingenix Suite Lab Data Management Software

