

## The paper-based QC process - 11 steps removed

