

## The paper-based QC process

