

Effective immediately, the LSCA Quality Manual has been replaced with the updated Agilent Quality Manual.

The URL link to the new manual is

[http://sharedoc.collaboration.agilent.com/sites/agilent-raqa/SD/Control of Quality Policy and Manual/Enterprise Docs/AQM01.pdf](http://sharedoc.collaboration.agilent.com/sites/agilent-raqa/SD/Control%20of%20Quality%20Policy%20and%20Manual/Enterprise%20Docs/AQM01.pdf)

Please update your web browser or any other documents to this new URL.

Questions? Submit them to enterprise.quality@agilent.com