



Agilent Technologies

August 31, 2009

If you need a copy of the Electronic Measurements Group Business Management System manual, please contact Ted Tucker, EMG Quality Manager, at ted_tucker@agilent.com.