

MANAGEMENT SYSTEM CERTIFICATE

Certificate no.:
CERT-08804-2006-AQ-HOU-ANAB

Initial certification date:
07 February, 1994

Valid:
08 August, 2024 – 07 August, 2027

This is to certify that the management system of
Micromeritics Instrument Corporation
4356 Communication Drive, Norcross, GA, 30093, USA

has been found to conform to the Quality Management System standard:
ISO 9001:2015

This certificate is valid for the following scope:
**Design, Manufacture and Service of Automatic and Manual Analytical Instruments which
Measure the Physical Characteristics of Powders and Solids**

Place and date:
Katy, TX, 06 August, 2024

For the issuing office:
DNV - Business Assurance
1400 Ravello Drive, Katy, TX, 77449-5164, USA



Sherif Mekkawy
Management Representative