

Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 9001:2015

This is to certify that: Automation Partnership
Grantham Close
York Way
Royston
SG8 5WY
United Kingdom

Holds Certificate Number: FS 533949

and operates a Quality Management System which complies with the requirements of ISO 9001:2015 for the following scope:

The design, development, manufacture and installation of automation systems for life-science industries, including the provision of servicing, technical support and consultation.



For and on behalf of BSI:

Andrew Launn - EMEA Systems Certification Operations and Compliance Director

Original Registration Date: 2008-07-03

Latest Revision Date: 2017-06-21

Effective Date: 2017-09-04

Expiry Date: 2020-09-03

Page: 1 of 1



003

...making excellence a habit.™