



MANAGEMENT SYSTEM CERTIFICATE

Certificate no.:
241476-2017-AQ-FIN-FINAS

Initial certification date:
12 June 2017

Valid:
28 October 2024 – 12 June 2026

This is to certify that the management system of

Sophion Bioscience A/S

Industriparken 39, 2750 Ballerup, Denmark

has been found to conform to the Quality Management System standard:

ISO 9001:2015

This certificate is valid for the following scope:

Development, Manufacturing, Sales and Customer Support of precision instruments for research and development within drug discovery.

Place and date:
Espoo, 28 October 2024

For the issuing office:
DNV - Business Assurance
Keilaranta 1, 02150 Espoo, Finland



Kimmo Haarala
Management Representative