



# MANAGEMENT SYSTEM CERTIFICATE

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

Initial certification date:  
12 June 2017

Valid:  
13 June 2026 – 12 June 2029

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, 04 June 2026



For the issuing office:  
DNV - Business Assurance  
Keilaranta 15B, 02150 Espoo, Finland

**Kimmo Haarala**  
Management Representative

Lack of fulfilment of conditions as set out in the Certification Agreement may render this Certificate invalid.

ACCREDITED UNIT: DNV Business Assurance Finland Oy Ab, Keilaranta 15B, 02150 Espoo, Finland - TEL: +358 10 292 4200. [www.dnv.fi/assurance](http://www.dnv.fi/assurance)