



CERTIFICATE



This is to certify that



InSphero AG

Wagistrasse 27a
8952 Schlieren
Switzerland

has implemented and maintains a **Quality Management System**.

Scope:

Development, production and distribution of new organotypic 3D Microtissues for toxicology and efficacy studies for chemistry, pharmaceuticals, cosmetics and nutritional applications Compound testing services using 3D Microtissues Development, production and distribution of laboratory consumables and instrument for 3D cell culture

Through an audit, documented in a report, it was verified that the management system fulfills the requirements of the following standard:

ISO 9001 : 2015

Certificate registration no.	495395 QM15
Valid from	2023-12-23
Valid until	2026-12-22
Date of certification	2023-11-03



DQS IS A MEMBER OF



DQS GmbH

Christian Gerling
Managing Director



Accredited Body: DQS GmbH, August-Schanz-Straße 21, 60433 Frankfurt am Main, Germany
The validity of this certificate can only be verified by the QR-code.

Certificate

DQS Holding has issued an IQNET recognized certificate that the organization:

InSphero AG
Wagistrasse 27a
CH-8952 Schlieren Switzerland

has implemented and maintains a
Quality Management System

for the following scope:

Development, production and distribution of new organotypic 3D Microtissues for toxicology and efficacy studies for chemistry, pharmaceuticals, cosmetics and nutritional applications Compound testing services using 3D Microtissues Development, production and distribution of laboratory consumables and instrument for 3D cell culture

which fulfils the requirements of the following Standard:

ISO 9001:2015

Issued on: **2023-12-23**
Expires on: **2026-12-22**

Registration Number: **DE-495395 QM15**




Alex Stoichitoiu
President of IQNet


Pedro Alves
Managing Director of IQNet

.....
This annex is only valid in connection with the original certificate number mentioned above.
.....

IQNET Members*:

AENOR Spain **AFNOR Certification** France **APCER** Portugal **CCC** Cyprus **CISQ** Italy **CQC** China **CQM** China **CQS** Czech Republic
Cro Cert Croatia **DQS Holding GmbH** Germany **EAGLE Certification Group** USA **FCAV** Brazil **FONDONORMA** Venezuela **ICONTEC**
Colombia **ICS** Bosnia and Herzegovina **INTECO** Costa Rica **IRAM** Argentina **JQA** Japan **KFQ** Korea **LSQA** Uruguay **MIRTEC** Greece
MSZT Hungary **Nemko AS** Norway **NSAI** Ireland **NYCE** México **PCBC** Poland **Quality Austria** Austria **SII** Israel **SIQ** Slovenia **SIRIM**
QAS International Malaysia **SQS** Switzerland **SRAC** Romania **TSE** Türkiye **YUQS** Serbia

* The list of IQNET Members is valid at the time of issue of this certificate. Updated information is available under www.iqnet-certification.com