



CERTIFICATO n. **316**
 CERTIFICATE No

SI CERTIFICA CHE L'ORGANIZZAZIONE
 WE HEREBY CERTIFY THAT THE ORGANIZATION

BIO-RAD LABORATORIES SRL

IT - 20090 SEGRATE (MI) - VIA CELLINI, 18/A

NELLE SEGUENTI UNITA' OPERATIVE / IN THE FOLLOWING OPERATIVE UNITS

IT - 20090 SEGRATE (MI) - VIA CELLINI 18/A

28109 ALCOBENDAS - MADRID EL SOTO DE LA MORALEJA - CALENDULA, 95 - MINIPARC II - EDIF. M

PT - 2610-156 AMADORA (LISBON) ALFRAGIDE ED. PRIME - AV. QUINTA GRANDE 53 - 3B -

GR - 11527 ATHENS AMPELOKIPI - 2-4 MESOGION AVE. (ATHENS TOWER)

HA ATTUATO E MANTIENE UN SISTEMA DI GESTIONE QUALITA' CHE E' CONFORME ALLA NORMA
 HAS IMPLEMENTED AND MAINTAINS A QUALITY MANAGEMENT SYSTEM WHICH COMPLIES WITH THE FOLLOWING STANDARD

UNI EN ISO 9001:2015

PER LE SEGUENTI ATTIVITA' / FOR THE FOLLOWING ACTIVITIES

SETTORE
 CODE

IAF 29, 19

Commercializzazione di strumenti e reagenti per la caratterizzazione di materiali biologici e sistemi
 per la ricerca e per applicazioni industriali.

Commercializzazione di sistemi, reagenti, materiali di controllo, dispositivi medici diagnostici in vitro e
 kits per la biologia molecolare, l' immunoematologia e la tipizzazione dei gruppi sanguigni.

Servizi di assistenza tecnica e specialistica al Cliente.

*Sales of instruments and reagents for biological material characterization and systems for research
 and industrial use.*

*Sale of systems, reagents, control materials, in-vitro diagnostics medical devices and kits for
 molecular biology, immunohematology and blood typing.*

Technical and specialized assistance to Customers.

IL PRESENTE CERTIFICATO E' SOGGETTO AL RISPETTO DEL REGOLAMENTO PER LA CERTIFICAZIONE DEI SISTEMI DI GESTIONE
 THE USE AND THE VALIDITY OF THE CERTIFICATE SHALL SATISFY THE REQUIREMENTS OF THE RULES FOR THE CERTIFICATION OF MANAGEMENT SYSTEMS

PRIMA EMISSIONE **23/06/1995**
 FIRST ISSUE
 DATA DELIBERA **27/03/2018**
 DECISION DATE
 DATA SCADENZA **08/04/2021**
 EXPIRY DATE
 EMISSIONE CORRENTE **19/09/2019**
 ISSUE DATE

CERTIQUALITY S.r.l. - IL PRESIDENTE
 Via G. Giardino 4 - 20123 MILANO (MI) - ITALY



SGQ n. 008 A SSI N. 007 G
 SGA n. 001 D SGE N. 001 M
 SCR n. 002 F ISP N. 006 O
 FSM n. 006 I EMAS N. 008 P
 PRD n. 008 B ITX N. 004 L
 DAP n. 003 H
 PRS n. 100 C

Membro degli Accordi di Mutuo
 Riconoscimento EA, IAF e ILAC
 Signatory of EA, IAF and ILAC Mutual
 Recognition Agreements

CISQ is a member of



IO Net, the association of the world's first class
 certification bodies, is the largest provider of management
 System Certification in the world.
 IO Net is composed of more than 30 bodies and counts
 over 150 subsidiaries all over the globe.

For information concerning the validity
 of the certificate, you can visit the site
www.certiquality.it

The validity this certificate depends
 on annual audit and on a complete
 review every three years of the
 Management System.



www.cisq.com

CISQ è la Federazione Italiana di Organismi di
 Certificazione dei sistemi di gestione aziendale.
 CISQ is the Italian Federation of management
 system Certification Bodies.



THE INTERNATIONAL CERTIFICATION NETWORK

CERTIFICATE

CISQ/CERTIQUALITY S.r.l.

has issued an IQNet recognised certificate that the organization:

BIO-RAD LABORATORIES SRL

IT - 20090 SEGRATE (MI) - VIA CELLINI, 18/A

for the following scope

Sales of instruments and reagents for biological material characterization and systems for research and industrial use.

Sale of systems, reagents, control materials, in-vitro diagnostics medical devices and kits for molecular biology, immunohematology and blood typing.

Technical and specialized assistance to Customers.

has implemented and maintains a

Quality Management System

which fulfills the requirements of the following standard

ISO 9001:2015

Issued on: **2018-03-27**

First issued on: **1995-06-23**

Expires on: **2021-04-08**

This attestation is directly linked to the IQNet Partner's original certificate and shall not be used as a stand-alone document

Registration number: **IT-1800**



Alex Stoichitoiu
President of IQNET



Ing. Claudio Provetti
President of CISQ

IQNet Partners*:

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 FCAV Brazil
FONDONORMA Venezuela ICONTEC Colombia Inspecta Sertifiointi Oy Finland INTECO Costa Rica
IRAM Argentina JQA Japan KFQ Korea MIRTEC Greece MSZT Hungary Nemko AS Norway NSAI Ireland
NYCE-SIGE México PCBC Poland Quality Austria Austria RR Russia SII Israel SIQ Slovenia
SIRIM QAS International Malaysia SQS Switzerland SRAC Romania TEST St Petersburg Russia TSE Turkey YUQS Serbia
IQNet is represented in the USA by: AFNOR Certification, CISQ, DQS Holding GmbH and NSAI Inc.