



BUREAU
VERITAS

PROMEGA ITALIA S.R.L

VIALE PIERO E ALBERTO PIRELLI, 6 - 20126 MILANO (344) - Italy

Certified site:

VIALE PIERO E ALBERTO PIRELLI, 6 - 20126 MILANO (344) - Italy

Bureau Veritas Certification Holding SAS - Italy Branch certifies that the Management System of the above organisation has been audited and found to be in accordance with the requirements of the management system standards detailed below

ISO 9001:2015

Scope of certification

Marketing of reagents and instruments for biomedical research and molecular diagnostics.

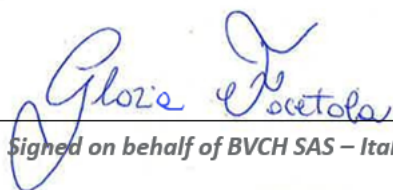
IAF : 29

Original Cycle Start Date:	19-March-2004
Expiry date of previous cycle:	29-June-2025
Certification / Recertification Audit date:	07-May-2025
Certification/Recertification Cycle Start date:	11-June-2025
Subject to the continued satisfactory operation of the organization's Management System, this certificate expires on:	29-June-2028

Certificate No. **IT343199**

Version: 1

Revision date: 11-June-2025


Signed on behalf of BVCH SAS – Italy Branch



02138

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

Certification body address: Viale Monza, 347 - 20126 Milano (MI) Italy
Local office: Viale Monza, 347 - 20126 Milano (MI) Italy

Further clarifications regarding the scope of this certificate and the applicability of the management system requirements may be obtained by consulting the organization.
To check the validity of this certificate please double click or scan QR CODE

