



BUREAU
VERITAS

PROMEGA ITALIA S.R.L.

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

Certified site:

Viale Piero e Alberto Pirelli, 6 - 20126 MILANO (MI) - Italy

Bureau Veritas Italia S.p.A. certifies that the Management System of the above organization has been audited and found to be in accordance with the requirements of the management system standards detailed below

ISO 14001:2015

Scope of certification

Sales of reagents and related equipment for biomedical research and molecular diagnostics.

IAF:29

Certificate issued in accordance with the Technical Regulation ACCREDIA RT-09

Original Cycle Start Date:	14-June-2024
Expiry date of previous cycle:	NA
Certification / Recertification Audit date:	13-May-2024
Certification/Recertification Cycle Start date:	14-June-2024
Subject to the continued satisfactory operation of the organization's Management System, this certificate expires on:	13-June-2027

Certificate No. **IT332931**

Version: 1

Revision date: 14-June-2024



 GLORIA FOCETOLA - Local Technical Manager



MS N°0009

Certification Body Address:
Bureau Veritas Italia S.p.A., Viale Monza, 347 - 20126 Milan, 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

