

**BUREAU VERITAS**  
Certification



## PROMEGA ITALIA SRL

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

**Certified site:**

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

*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 9001:2015

*Scope of certification*

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

IAF sector(s): **29**

Original cycle start date: **19 March 2004**

Expiry date of previous cycle: **NA**

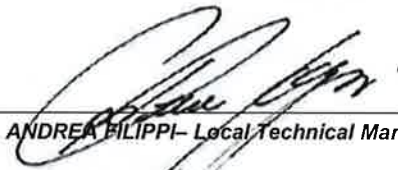
Certification / Recertification Audit date: **04 June 2019**

Certification / Recertification cycle start date: **14 June 2019**

Subject to the continued satisfactory operation of the organization's Management System, this certificate expires on: **29 June 2022**

Certificate No. - Version: **216909- 1**

Revision date: **14 June 2019**

  
**ANDREA FILIPPI - Local Technical Manager**

Certification body address:  
Bureau Veritas Italia S.p.A., Viale Monza, 347 - 20126 Milano, Italia

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 this certificate validity please refer to the website [www.bureauveritas.it](http://www.bureauveritas.it)



SGQ N° 009A

Member degli Accordi di Mutuo Riconoscimento EA, IAF e ILAC  
Signatory of EA, IAF and ILAC mutual Recognition Agreements

