

BUREAU VERITAS
Certification



PROMEGA ITALIA SRL

Registered and Operative Site:

Via Decembrio Pier Candido, 28 – 20137 MILANO (MI) - ITALY

Bureau Veritas Italia spa certify 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

Standard

ISO 9001:2015

Scope of certification

Distribution of reagents and related equipment for biomedical research.

EA Sector(s): **29**

Certification cycle start date: **09 June 2016**

Subject to the continued satisfactory operation of the organization's Management System, this certificate expires on: **29 June 2019**

Original certification date: **19 March 2004**

Certificate No. **216909**

Version N. **1**

Revision date: **09 June 2016**


ANDREA FILIPPI – Local Technical Manager

Certification body address:

Bureau Veritas Italia spa, Via Miramare, 15, 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



SGQ	N° 009A	PRS	N° 076C
SGA	N° 008D	SGE	N° 009M
PRD	N° 009B	EMAS	N° 004P
SCR	N° 008F	GHG	N° 008D
FSMS	N° 003T	ISP	N° 006E

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