

BUREAU VERITAS
Certification



Certification

Awarded to

PROMEGA CORPORATION (AUSTRALIA)

SUITE W3A, LEVEL 3, BUILDING 2, SYDNEY CORPORATE PARK, 75 O'RIORDAN
STREET, ALEXANDRIA, NSW, 2015

SUITE 4.05, 737 BURWOOD ROAD, HAWTHORN, VIC, 3123

AUSTRALIA

Bureau Veritas 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 indicated below

STANDARD

ISO 9001:2015

SCOPE OF SUPPLY

ALEXANDRIA, NSW: IMPORTATION, STOCKHOLDING, SALES AND TECHNICAL SUPPORT OF BIOLOGICAL RESEARCH PRODUCTS. IMPORTATION AND MANAGEMENT OF CONTRACT SERVICING OF RELATED EQUIPMENT

HAWTHORN, VIC: SALES AND TECHNICAL SUPPORT

Original Approval Date: **28 January 1998**

Subject to the continued satisfactory operation of the organisation's Management System,

this certificate is valid until: **31 January 2022**

To check the validity of this certificate please call tel. **1800 855 190**

Further clarification regarding the scope of this certificate and the applicability of the Management System requirements may be obtained by consulting the organisation.

Certificate Number: **AU003405-1**

Date: **22 January 2019**

Andrew Mortimore
Vice President – I&F Pacific Region

Managing office: Bureau Veritas Pty Ltd, 3/435 Williamstown Road,
Port Melbourne, Victoria, 3207

Issuing office: Bureau Veritas Pty Ltd, 3/435 Williamstown Road,
Port Melbourne, Victoria, 3207

JAS-ANZ



www.jas-anz.org/register



**BUREAU
VERITAS**