

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
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

**IMPORTATION, STOCKHOLDING, SALES AND TECHNICAL SUPPORT OF
BIOLOGICAL RESEARCH PRODUCTS. IMPORTATION AND MANAGEMENT OF
CONTRACT SERVICING OF RELATED EQUIPMENT. SERVICING OF MAXWELL AND
GLOMAX INSTRUMENTS**

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 2028**

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: **AU006350-1**

Date: **10 January 2025**

Andrew Mortimore
VP/CE CIF Pacific Region

Managing office: Bureau Veritas Pty Ltd, Level 11, 500 Collins Street,
Melbourne, Victoria, 3000, Australia

Issuing office: Bureau Veritas Pty Ltd, Level 11, 500 Collins Street,
Melbourne, Victoria, 3000, Australia

JAS-ANZ



www.jas-anz.org/register



**BUREAU
VERITAS**