

**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 2025**

*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:* **AU004867-1**

*Date:* **24 January 2022**

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](http://www.jas-anz.org/register)



**BUREAU  
VERITAS**