



BUREAU
VERITAS

Bureau Veritas Certification

Certification

Awarded to

PROMEGA CORPORATION

2800 WOODS HOLLOW ROAD, MADISON, WI, 53711, United States

This is a multisite certificate, additional site(s) are listed on the next page(s)

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 18385:2016

Scope of Supply

THE PRODUCTION OF PRODUCTS USED IN THE COLLECTION, STORAGE, AND ANALYSIS OF BIOLOGICAL MATERIAL FOR FORENSIC DNA PURPOSES, AND THE REQUIREMENT FOR MANUFACTURERS TO MINIMIZE THE RISK OF OCCURRENCE OF DETECTABLE HUMAN NUCLEAR DNA CONTAMINATION IN PRODUCTS USED BY THE GLOBAL FORENSIC COMMUNITY

Original approval date: 03-05-2017

Subject to the continued satisfactory operation of the organisation's Management System, this certificate is valid until: 02-05-2029

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 No.: **AU006948**

Version: 1

Date: **20-04-2026**

Andrew Mortimore
VP/CE CIF Pacific Region

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

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





BUREAU
VERITAS

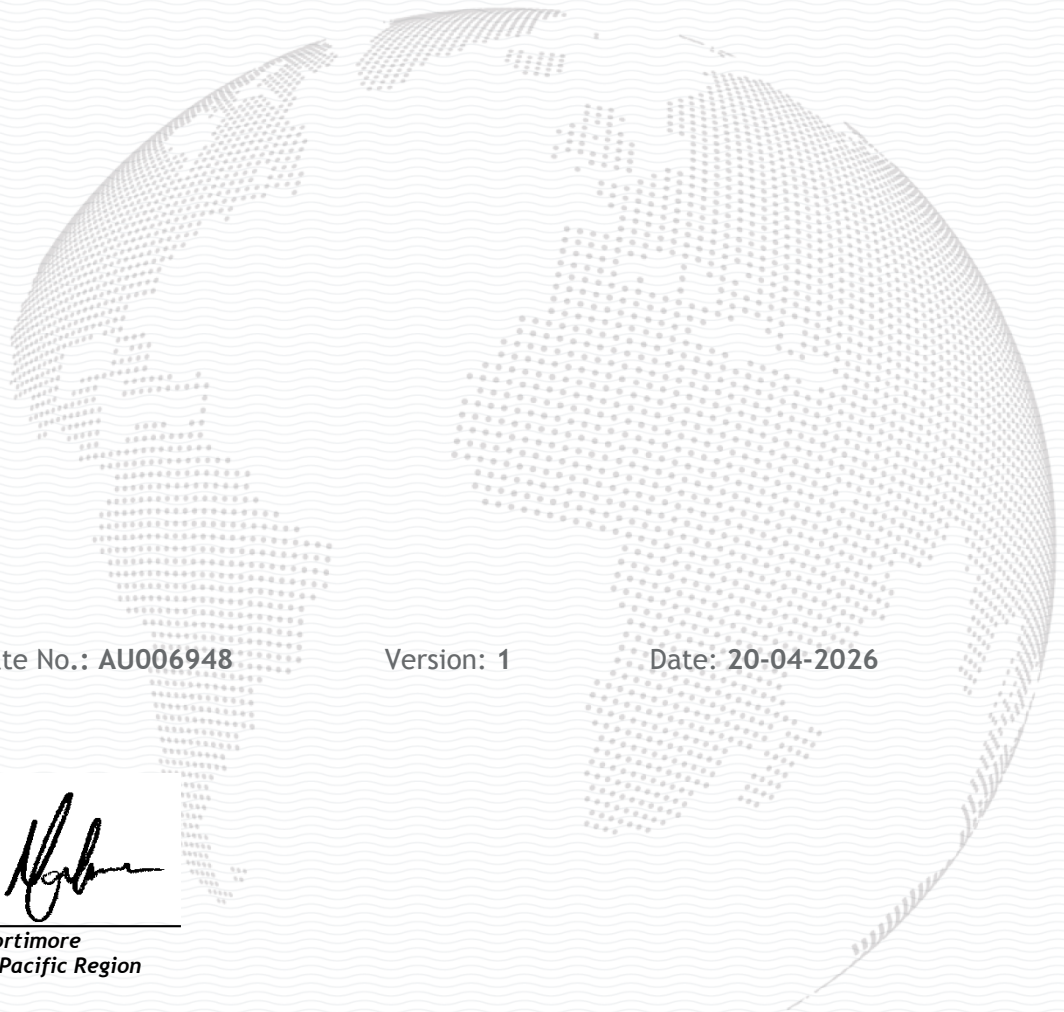
PROMEGA CORPORATION

STANDARD

ISO 18385:2016

Scope of supply

Site Name	Site Address	Site Scope
Head Office	2800 WOODS HOLLOW ROAD, MADISON, WI, 53711, United States	THE PRODUCTION OF PRODUCTS USED IN THE COLLECTION, STORAGE, AND ANALYSIS OF BIOLOGICAL MATERIAL FOR FORENSIC DNA PURPOSES, AND THE REQUIREMENT FOR MANUFACTURERS TO MINIMIZE THE RISK OF OCCURRENCE OF DETECTABLE HUMAN NUCLEAR DNA CONTAMINATION IN PRODUCTS USED BY THE GLOBAL FORENSIC COMMUNITY
BioPharmaceutical Technology Center	5445 East Cheryl Parkway, Fitchburg, Wisconsin, 53711, United States	
Kepler Center	5455 Nobel Drive, Fitchburg, Wisconsin, 53711, United States	



Certificate No.: **AU006948**

Version: **1**

Date: **20-04-2026**

Andrew Mortimore
VP/CE CIF Pacific Region

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

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

