

BUREAU VERITAS
Certification



Certification

Awarded to

PROMEGA CORPORATION

BIOPHARMACEUTICAL TECHNOLOGY CENTER-

5445 E. CHERYL PARKWAY, FITCHBURG WI 53711

FEYNMAN CENTER

2780 WOODS HOLLOW ROAD, FITCHBURG WI 53711

KEPLER CENTER

5455 NOBEL DRIVE, FITCHBURG, WI 53711

UNITED STATES OF AMERICA

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: **3 May 2017**

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

this certificate is valid until: **02 May 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 Number: **AU006498-1**

Date: **12 March 2026**

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



**BUREAU
VERITAS**