

BUREAU VERITAS
Certification



Certification

Awarded to

PROMEGA CORPORATION

**2800 WOODS HOLLOW ROAD, MADISON, 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, INCLUDING REAGENTS AND CONSUMABLES,
USED BY THE GLOBAL FORENSIC COMMUNITY FOR THE PREPARATION AND
ANALYSIS OF BIOLOGICAL MATERIAL USED FOR FORENSIC PURPOSES, WHILE
MINIMIZING THE RISK OF DETECTABLE HUMAN DNA CONTAMINATION**

Original Approval Date: **3 March 2017**

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

this certificate is valid until: **02 May 2023**

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: **AU003930-1**

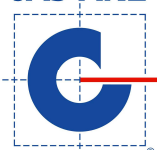
Date: **24 March 2020**

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**