



MANAGEMENT SYSTEM CERTIFICATE

Certificate no.:
178611-2015-AQ-VNM-JAS-ANZ

Initial certification date:
16 July 2009

Valid:
17 July 2024 – 16 July 2027

This is to certify that the management system of

PMAS CO., LTD.

Lot A-2A-CN, My Phuoc 3rd Industrial Park, My Phuoc Ward, Ben Cat Town, Binh Duong Province, Vietnam

has been found to conform to the Quality Management System standard:

ISO 9001:2015

This certificate is valid for the following scope:

Precision Machining Components and Assembly Services.

Place and date:
Singapore, 09 May 2024

For the issuing office:
DNV - Business Assurance
16 Science Park Drive, DNV Technology
Centre, Singapore 118227, Singapore



Siang Hooi Silas Ding
Management Representative

Lack of fulfilment of conditions as set out in the Certification Agreement may render this Certificate invalid.

Accreditation by the joint accreditation system of Australia and New Zealand (www.jas-anz.org/register)
ACCREDITED UNIT: DNV Business Assurance Australia Pty Limited, Level 7, 124 Walker Street, North Sydney, NSW, 2060, Australia
TEL: +61 2 9900 9500. <https://www.dnv.com.au/assurance/>