

Certificate IATF 0511849

Certificate MY17/02441

The management system of

# Paramit Malaysia Sdn. Bhd.

IATF USI: TA4E4L

PMT 1199 (Plot 372), Penang Science Park, Lorong Perindustrian Bukit Minyak 21, Taman Perindustrian Bukit Minyak, Simpang Ampat, 14100 Seberang Perai Tengah, Pulau Pinang, Malaysia

has been assessed and certified as meeting the requirements of

## IATF 16949:2016

Edition 1

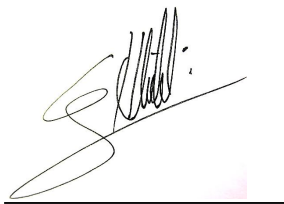
For the following Scope

Manufacture of Electronic Modules and Electromechanical Products

### EXCLUSIONS : 8.3 Product Design

36 Month certification is valid from 17 April 2024 until 16 April 2027 and remains valid subject to satisfactory surveillance audits.

Version no. 4. Current version updated 17 April 2024



Authorised by

Gatot Hartanto

Veto Power Authority

Contracted Office : SGS United Kingdom Ltd, Station Road, Oldbury, B69 4LN, UK.

Email : [Neil.Hall@sgs.com](mailto:Neil.Hall@sgs.com)



This document is an authentic electronic certificate for Client' business purposes use only. Printed version of the electronic certificate are permitted and will be considered as a copy. This document is issued by the Company subject to SGS General Conditions of certification services available on [Terms and Conditions | SGS](#). Attention is drawn to the limitation of liability, indemnification and jurisdictional clauses contained therein. This document is copyright protected and any unauthorized alteration, forgery or falsification of the content or appearance of this document is unlawful.

