

Certificate MY96/07435

The management system of

# Paramit Malaysia Sdn. Bhd.

# SGS

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  
**ISO 9001:2015**

For the following activities  
Manufacture of Electronic Modules and Electromechanical Products

This certificate is valid from 17 April 2024 until 16 April 2027 and remains valid subject to satisfactory surveillance audits.

Issue 16. Certified since 20 June 1996



Authorised by  
Jonathan Hall  
Global Head - Certification  
Services

SGS United Kingdom Ltd  
Rossmore Business Park, Ellesmere Port, Cheshire, CH65 3EN, UK  
t +44 (0)151 350-6666 - [www.sgs.com](http://www.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.

