



Certificate MY04/61076

The management system of

Paramit Malaysia Sdn. Bhd.

PMT 1199 (Plot 372), Penang Science Park
Lorong Perindustrian Bukit Minyak 21, Taman Perindustrian Bukit Minyak
14100 Simpang Ampat, Seberang Prai Tengah, Penang
MALAYSIA

has been assessed and certified as meeting the requirements of

ISO 14001:2015

For the following activities

**Contract Manufacturing Service Provider of Printed Circuit Assemblies
and System Box Builds.**

This certificate is valid from 20 January 2022 until 19 January 2025 and
remains valid subject to satisfactory surveillance audits.
Recertification audit due a minimum of 60 days before the expiration date.
Issue 11. Certified since 20 January 2004



Authorised by



0005

SGS United Kingdom Ltd
Rossmore Business Park, Ellesmere Port, Cheshire, CH65 3EN, UK
t +44 (0)151 350-6666 f +44 (0)151 350-6600 www.sgs.com

21HC 14001:2015_0421

Page 1 of 1



This document is a Web version of SGS certificate for electronic use exclusively. It shall only be available by clicking on SGS Certification Mark which has been posted on Your website. It shall not be printed in anyway. This document is copyright protected. No content or appearance may be reproduced without the express written permission of SGS. Any misuse, alteration, forgery or falsification is unlawful.