

Certificate MY12/00914

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 Perai Tengah, Pulau Pinang, Malaysia

has been assessed and certified as meeting the requirements of

ISO 13485:2016

EN ISO 13485:2016

For the following activities

Contract Manufacture of Printed Circuit Assemblies, System Box Builds for Active Medical Device (Non-implantable) and Finished Diagnostic & Instrumentation Devices in accordance with Customer Specified Requirements.

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

Issue 7. Certified since 17 April 2012

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



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.

