

Certificate MY04/61076

The management system of

# Tecan CDMO Solutions PN 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 2025 until 20 January 2028 and remains valid subject to satisfactory surveillance audits.

Issue 13. Certified since 20 January 2004

*L. Moran*

Authorised by

Liz Moran  
Business Manager

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.

