

Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 13485:2016 & EN ISO 13485:2016

This is to certify that:

KANEKA EUROGENTEC S.A.
Rue Bois Saint-Jean, 5
Liège Science Park
Seraing
B-4102
Belgium

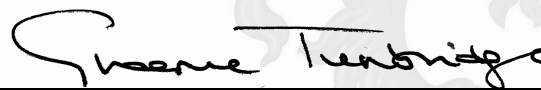
Holds Certificate Number:

MD 638600

and operates a Quality Management System which complies with the requirements of ISO 13485:2016 & EN ISO 13485:2016 for the following scope:

Contract manufacturing of synthetic oligonucleotides in support of in vitro diagnostic and related applications.
Production d'oligonucléotides synthétiques en support au diagnostic in vitro et aux applications associées.

For and on behalf of BSI:



Graeme Tunbridge, Senior Vice President Medical Devices

Original Registration Date: 2015-07-23

Latest Revision Date: 2023-11-21

Effective Date: 2023-12-23

Expiry Date: 2026-12-22



Page: 1 of 1

...making excellence a habit.™