

Certificate of Registration

QUALITY MANAGEMENT SYSTEM - ISO 9001:2015

This is to certify that:

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

Holds Certificate Number:

FS 638601

and operates a Quality Management System which complies with the requirements of ISO 9001:2015 for the following scope:

Development, production and sales of products and services for research and development in life sciences.

Développement, production et vente de produits et services pour la recherche et le développement dans les sciences de la vie.

For and on behalf of BSI:

Matt Page, Managing Director Assurance - UK & Ireland

Original Registration Date: 2015-07-27

Effective Date: 2023-12-23

Latest Revision Date: 2023-11-15

Expiry Date: 2026-12-22



Page: 1 of 2

...making excellence a habit.™

Certificate No: FS 638601

Location

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

KANEKA EUROGENTEC S.A.
UK Branch
Unit 4, Albany Court
Albany Park
Camberly
GU16 7QR
United Kingdom

Registered Activities

Development, production and sales of products and services for research and development in life sciences.

Développement, production et vente de produits et services pour la recherche et le développement dans les sciences de la vie.

Production and Quality Control of products for research and development in Life sciences.

Production et Contrôle Qualité de produits pour la recherche et développement dans les sciences de la vie.



Original Registration Date: 2015-07-27

Latest Revision Date: 2023-11-15

Effective Date: 2023-12-23

Expiry Date: 2026-12-22

Page: 2 of 2

This certificate was issued electronically and remains the property of BSI and is bound by the conditions of contract.

An electronic certificate can be authenticated [online](#).

Printed copies can be validated at www.bsigroup.com/ClientDirectory

Information and Contact: BSI, Kitemark Court, Davy Avenue, Knowlhill, Milton Keynes MK5 8PP. Tel: + 44 345 080 9000
BSI Assurance UK Limited, registered in England under number 7805321 at 389 Chiswick High Road, London W4 4AL, UK.
A Member of the BSI Group of Companies.