

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:

Andrew Launn, EMEA Systems Certification Director

Original Registration Date: 2015-07-27

Effective Date: 2020-12-23

Latest Revision Date: 2020-12-10

Expiry Date: 2023-12-22



003

Page: 1 of 2

...making excellence a habit.™

Certificate No: FS 638601

Location	Registered Activities
KANEKA EUROGENTEC S.A. Rue Bois Saint-Jean, 5 Liège Science Park Seraing B-4102 Belgium	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.
KANEKA EUROGENTEC S.A. UK Branch Unit 4, Albany Court Albany Park Camberly GU16 7QR United Kingdom	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.

Original Registration Date: 2015-07-27

Latest Revision Date: 2020-12-10

Effective Date: 2020-12-23

Expiry Date: 2023-12-22

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.