

CERTIFICAT
CERTIFICATE OF REGISTRATION
N° 15217 - Revision 1

Le LNE certifie que le système de management de la qualité développé par
LNE certifies that the quality management system developed by

EUROGENTEC S.A.
LIEGE science park, Rue Bois Saint-Jean, 5
4102 SERAING BELGIUM

pour les activités
for the activities

Production et vente d'oligonucléotides de diagnostic in vitro

Production and sales of in vitro diagnostic oligonucleotides

réalisées sur le(s) site(s) de
performed on the location(s) of

EUROGENTEC S.A.
LIEGE Science Park - Rue Bois Saint Jean, 5 - 4102 SERAING - BELGIUM

est conforme aux exigences des normes internationales
complies with the requirements of the international standards

ISO 13485 : 2003

Début de validité / Effective date : December 23rd, 2011 (included)
Valable jusqu'au / Expiry date : December 22nd, 2014 (included)
Etabli le / Issued on : November 28th, 2011



For the General Director
Laurence DAGALLIER
Deputy Director