

Certificate

Standard **ISO 9001:2015**

Certificate Registr. No. 0.04.08238

Certificate Holder: **PROMEGA BIOTECH IBERICA, S.L.**
Avda. de Bruselas, 5-3ª planta
28108 Alcobendas (Madrid)
Spain

Scope: The distribution, sales and after-sales service of products and equipment for detecting and extracting nucleic acids and proteins.

Proof has been furnished by means of an audit that the requirements of ISO 9001:2015 are met.

Validity: The certificate is valid from 2024-02-10 until 2027-02-09.
First certification 2009

2024-01-10



TÜV Rheinland Ibérica Inspection,
Certification & Testing S.A.
Garrotxa, 10-12 – E-08820 El Prat de
Llobregat

