

Certificate

Standard **ISO 14001:2015**

Certificate Registr. No. 3.00.21034

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 14001:2015 are met.

Validity: The certificate is valid from 2022-01-31 until 2025-01-30.
First certification 2022

2022-03-09



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