

# 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 2025-01-31 until 2028-01-30.  
First certification 2022.

2025-01-21

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



**TÜVRheinland®**

www.tuv.com