



BUREAU
VERITAS

Bureau Veritas Certification

Certification

Awarded to

PID ENG & TECH, S.L.

**CL FRANCISCO GERVÁS 11 POL IND ALCOBENDAS - 28108 - ALCOBENDAS - MADRID
- SPAIN**

Bureau Veritas certification certifies that the Management System has been audited and found to be in accordance with the requirements of standard:

STANDARD

ISO 14001:2015

Scope of certification:

**DESIGN, ASSEMBLY, OPTIMIZATION AND TRADING OF SCIENTIFIC EQUIPMENT AND
BENCH SCALE PILOT PLANTS FOR PHYSICAL AND CHEMICAL PROCESSES.**

Certificate Number:	ES145933 - 1
Original approval date:	07-01-2007
Certification/Renovation Audit:	24-05-2024
Expiry date of previous cycle:	25-07-2024
Effective date:	26-07-2024
Certificate expiration date:	25-07-2027

This certificate is valid, subject to the general and specific terms and conditions of certification services

Bureau Veritas Iberia S.L.
C/ Valportillo Primera 22-24, 28108 Alcobendas - Madrid, España

