

  
**RI.  
SE**  
**CERTIFIKAT****ISO 14001**

Härmed intygas att/This is to certify that

**Promega Biotech AB**

Finnboda Varvsväg 16, 131 73 NACKA, SWEDEN

har ett miljöledningssystem som uppfyller kraven enligt SS-EN ISO 14001:2015 vad gäller  
has an environmental management system that fulfils the requirements of SS-EN ISO 14001:2015 with respect to

Försäljning och marknadsföring av produkter  
för forskning och utveckling inom Life Science

Sales and marketing of products for research  
and development in Life Science

Ursprungligen utfärdat/Originally issued	2019-12-10
Giltigt till och med/Expiry date	2022-12-09
Beslutsdatum/Decision date	2019-12-10

Dag Sjöholm

**RISE Research Institutes of Sweden AB** | Certification  
Box 857, SE-501 15 Borås, Sweden  
☎ +46 10-516 50 00  
certifiering@ri.se | www.ri.se



Certifikat/Certificate C000962 | version/issue 1 | 2019-12-10

This document is the property of RISE and may not be reproduced other than in full, except with the prior written approval by RISE.



Sida/Page 1(1)