



CERTIFIKAT

ISO 9001

Härmed intygas att/This is to certify that

Promega Biotech AB

Finnboda Varvsväg 16, 131 73 NACKA, SWEDEN

har ett kvalitetsledningssystem som uppfyller kraven enligt SS-EN ISO 9001:2015 vad gäller
has a quality management system that fulfils the requirements of SS-EN ISO 9001:2015 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	2010-12-20
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 5000
certifiering@ri.se | www.ri.se



Certifikat/Certificate 5340 | version/issue 6 | 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.

