



Management Service

CERTIFICATE

Certificate Registration No.: **12 100 54414 TMS** / Order No.: **707067170**

**The Certification Body
of TÜV SÜD Management Service GmbH**

certifies that the organization



Reaction Biology Europe GmbH
Engesserstr. 4
79108 Freiburg im Breisgau
Germany

for the scope

**Development, production and sale of
recombinant protein products, biochemical,
cellular and immuno-oncological assay services,
and in vivo testing services for the
development of new drugs to treat diseases**

has established and applies a Quality Management System.

An audit was performed and has furnished proof
that the requirements according to

DIN EN ISO 9001:2015

are fulfilled.

The certificate is valid from **2026-02-22** until **2029-02-21**.

Fred Wenke
Head of Certification Body
Munich, 2026-02-13



CERTIFICAT

CERTIFICADO

СЕРТИФИКАТ

認證證書

CERTIFICATE

ZERTIFIKAT