



Management Service

CERTIFICATE

The Certification Body
of TÜV SÜD Management Service GmbH

certifies that



Reaction Biology Europe GmbH

Engesserstr. 4
79108 Freiburg im Breisgau
Germany

has established and applies
a Quality Management System for

**Development, production and sale of
recombinant protein products, biochemical and
cellular assay services as well as in vivo testing services for
the development of new drugs targeting diseases,
for example in oncology.**

An audit was performed, Order No. **707067170**.

Proof has been furnished that the requirements
according to

ISO 9001:2015

are fulfilled.

The certificate is valid from **2020-11-26** until **2023-08-09**.

Previous certificate valid until 2020-08-09.

Certificate Registration No.: **12 100 54414 TMS**.

Product Compliance Management
Munich, 2020-11-27



CERTIFICAT

CERTIFICADO

СЕРТИФИКАТ

認證證書

CERTIFICATE

ZERTIFIKAT