



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