

CERTIFICATE

The certification body IFTA AG certifies

Jena Bioscience GmbH
Löbstedter Straße 71
07749 Jena
Germany

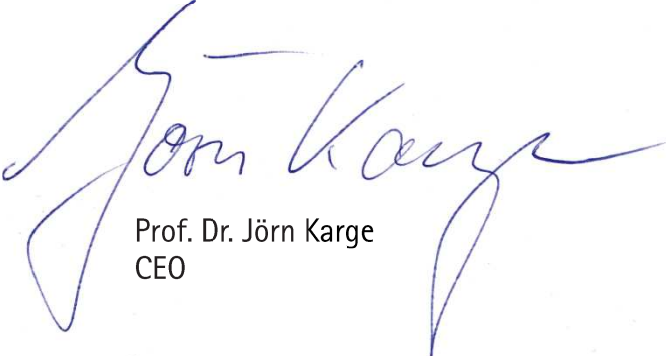
the conformity of the introduced quality management system for the field of development, production and distribution of reagents for life-sciences with the standard

DIN EN ISO 9001:2015

Start of validity:	13.03.2020
End of validity:	12.03.2023
Report and certificate number:	ICV03597 034 19
The certificate consists of	2 pages

This certificate includes an annual examination of the QMS by IFTA AG according to the specified standard.

Berlin, 02.03.2020



Prof. Dr. Jörn Karge
CEO

