

CERTIFICATE

The certification body IFTA AG certifies

Jena Bioscience GmbH
Löbstedter Straße 71
07749 Jena
Germany

the conformity of the introduced environmental management system with the standard

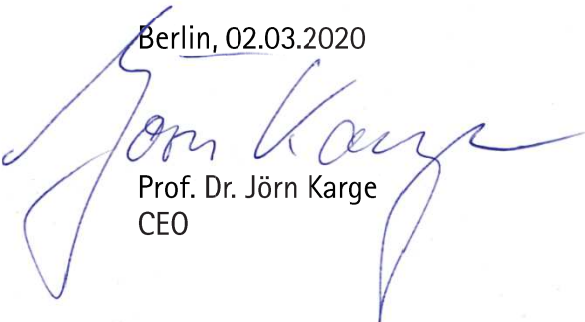
DIN EN ISO 14001:2015

for the environmentally relevant processes in the entire field of responsibility of the company, which relate to development, production and distribution of reagents for life-sciences.

Start of validity:	13.03.2020
End of validity:	12.03.2023
Report and certificate number:	ICV03597 534 19
The certificate consists of	2 pages

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

Berlin, 02.03.2020



Prof. Dr. Jörn Karge
CEO

