

# CERTIFICATE

This is a sub certificate and only valid  
in connection with the main certificate

ICC3597 034 11

The certification body confirms to

**Jena Bioscience GmbH**  
**Löbstedter Straße 80**  
**07749 Jena**  
**Germany**

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

**DIN EN ISO 9001:2008-12**

Start of validity:	01.12.2011
End of validity:	30.11.2014
Report and certificate number:	IC03214
The certificate consists of	1 page

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

Berlin, 01.12.2011

gez. Dr. Jörn Karge  
CEO



Deutscher  
Akkreditierungs  
Rat  
DGA-ZM-4453.00