



**Orion Registrar, Inc.**  
**Thorough and Fair Auditing**

## Certificate of Certification

*This is to certify the Quality Management System of:*

**LI-COR, Inc.**  
**4647 Superior Street**  
**Lincoln, NE 68504 USA**

*Has been assessed by Orion Registrar and found to be in compliance with the following Quality Standard:*

**ISO 9001:2015**

*The Quality Management System is applicable to:*

**Design, Manufacture, Market and Service of  
Biotechnology and Environmental Scientific  
Instruments and Reagents**

*The Certification period is from*

**June 1, 2017 to September 4, 2018**

*This certification is subject to the company maintaining its system to the required standard, and applicable exceptions, which will be monitored by Orion.*

Client ID: 430

Certificate ID: 1009085



  
*Paul M. Burck, President*      *June 1, 2017*  
Date

