



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 May 31, 2020

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: 1011761




Paul M. Burck, President *October 9, 2017*
Date

