

# CERTIFICATE of CALIBRATION for LI-COR SENSOR

Quantum Sensor  
Model Number: LI-190

Serial Number: **Q45044**

Calibration Date: **March 9, 2011**  
Manufacture Date: **February 9, 2011**

## Calibration Constant(s):

---

Output: 7.13 microamps per 1000  $\text{umol s}^{-1} \text{m}^{-2}$

## For use with LI-COR handheld meters and loggers:

---

Multiplier:  $-140.28 \text{ umol s}^{-1} \text{m}^{-2}$  per microamp

## If this is an LI-190-BL (3-wire bare leads):

---

Multiplier:  $140.28 \text{ umol s}^{-1} \text{m}^{-2}$  per microamp

## For use with LI-COR 2290 (604 ohm) millivolt Adapter:

---

Multiplier:  $-232.25 \text{ umol s}^{-1} \text{m}^{-2}$  per millivolt

## If this is an SL or SMV sensor:

---

Multiplier:  $-200 \text{ umol s}^{-1} \text{m}^{-2}$  per millivolt

**IMPORTANT:** Read the appropriate instruction manual ([www.licor.com/Quantum-manual](http://www.licor.com/Quantum-manual)) before using this sensor

**IMPORTANT:** It is recommended that sensors be recalibrated every two years after field deployment.

Calibration Technician: John Brooks

Calibration standard used: Working standard lamp 811C

Calibration traceable to the National Institute of Standards and Technologies (NIST) through NIST-calibration lamp number F-226.  
(NBS Test No. 534/237689-86)

**LI-COR**<sup>®</sup>

LI-COR Biosciences, Global Headquarters • 4647 Superior Street • Lincoln, NE 68504 USA  
Phone: +1-402-467-3576 • Fax: +1-402-467-2819 • Toll-free: 800-447-3576 (USA & Canada)  
[envsales@licor.com](mailto:envsales@licor.com) • [envsupport@licor.com](mailto:envsupport@licor.com) • [www.licor.com](http://www.licor.com)  
Manual: [www.licor.com/Quantum-manual](http://www.licor.com/Quantum-manual)