

Data Acquisition Module and Data Retention Module Specifications

The Data Acquisition Module connects digital, analog, and SDI-12 sensors to a LI-COR eddy covariance system. The Data Retention Module provides data logging and smart system controls that prioritize sensors if the power supply is compromised. This document describes the specifications of the two modules.

Data Acquisition Module (DAqM)

Configuration interface: Windows® and macOS graphical programming

Analog inputs:

Single-ended inputs: Up to 18 or

Differential inputs: Up to 9 (each differential input occupies two single-ended inputs)

With built-in excitation voltage: 2, 5 V, 25 mA

With selectable current or voltage: 2

Resolution: 24 bit

SDI-12 inputs

Digital inputs/outputs: 4

Output: 5 V logic

Input: 5 V or 3.3 V logic

Switched voltage outputs: 3

Output voltage: 9 to 30 VDC

Output current: Up to 1 A, with current sensing

12V excitation voltage output: 1

Up to 200 mA

Sampling rate: 1 measurement per second

Communications protocol: RS-485 input and output

Power requirements: +9 to 32 VDC

Power consumption: 0.7 to 1 W

Dimensions: 17 cm × 10.9 cm × 7 cm (W × H × D)

Weight: 0.52 kg

Data Retention Module (DRM)

Communication protocol: RS-485 input and output

Configurable power outputs: 4

Unregulated power output: 1 (power out equal to power supplied; 9 to 32 VDC)

Regulated power output: 3 (24 VDC maximum)

Power requirements: +9 to 32 VDC

Data storage: 32 MB; Over one month of backup biomet data collection, typically.

Dimensions: 17 cm × 10.9 cm × 7 cm (W × H × D; not including battery)

Weight: 0.52 kg (not including battery)

Backup Battery:

Type: Lead Acid

Output: 12 VDC; 7.2 Amp Hour



LI-COR Biosciences

4647 Superior Street
Lincoln, Nebraska 68504
Phone: +1-402-467-3576
Toll free: 800-447-3576 (U.S. and
Canada)
envsales@licor.com

LI-COR Distributor Network:
www.licor.com/env/distributors

©2018 LI-COR, Inc. All rights reserved.
982-17393 03/18

Regional Offices

LI-COR Biosciences GmbH
Siemensstraße 25A
61352 Bad Homburg
Germany
Phone: +49 (0) 6172 17 17 771
envsales-gmbh@licor.com

LI-COR Biosciences UK Ltd.
St. John's Innovation Centre
Cowley Road
Cambridge
CB4 0WS
United Kingdom
Phone: +44 (0) 1223 422102
envsales-UK@licor.com

Specifications subject to change.
LI-COR is an ISO 9001 registered
company. LI-COR is a registered
trademark of LI-COR, Inc. in the
United States and other countries.
All other trademarks belong to their
respective owners. For patent
information, visit
www.licor.com/patents.