

LI-7000 and LI-7500 Comparison

Specifications* for the LI-7000 Closed Path and LI-7500 Open Path CO₂/H₂O Analyzers



LI-7000



LI-7500

Gas Sample Path	Closed	Open
Analyzer Type	Differential NDIR	Absolute NDIR
Detector	Lead Selenide	Lead Selenide
Optical Bench	0.95 cm D × 15.2 cm L, 10.86 cm ³	1 cm D × 12.5 cm L optical path
Operating Temperature	0 to 50 °C	-25 to 50 °C (-40 to 50 °C optional range verified on request)
Power Consumption	15-40 W	10W in steady state, 30W during warm up
Pump	Yes	Not Required
Max. Cell Pressure	Capable of 10 liters/min. with minimal pressure drop	N/A
Pressure Sensor	Yes	Yes
Heat Exchangers	Yes	N/A
Tubing	External only	None
Weatherproof	No	Yes
Cleanable Optics	Yes	Yes
Windows® Interface	Yes	Yes
DACs	Four, updated at 600 Hz	Two, updated at 300 Hz
Digital Output	RS-232, USB up to 50 Hz	RS-232 to 20 Hz; SDM to 40 Hz
Bandwidth	0.025 to 20 Hz	5, 10 or 20 Hz
CO₂ Range	0-3,000 ppm	0-3,000 ppm
Zero Drift - max	±0.3 ppm/°C	±0.3 ppm/°C
Span Drift - max	±0.2% of reading/°C	±0.1% of reading/°C
Signal Noise	78 ppb at 5 Hz typical	80 ppb at 5 Hz typical
H₂O Range	0-60 mmol/mol	0-60 mmol/mol
Zero Drift - max	±0.02 mmol/mol/°C	±0.05 mmol/mol/°C
Span Drift - max	±0.4% of reading/°C	±0.3% of reading/°C
Signal Noise	0.004 mmol/mol at 5 Hz typical	0.003 mmol/mol at 5 Hz typical

*All specifications are subject to change without prior notice.

LI-COR®

Biosciences

LI-COR, Inc. • 4421 Superior Street • P.O. Box 4425 • Lincoln, NE 68504. U.S.A.
 Toll Free: 1-800-447-3576 (U.S. & Canada) • Phone: 402-467-3576 • FAX: 402-467-2819
 In Germany - LI-COR GmbH: +49 (0) 6172 17 17 771
 E-mail: envsales@licor.com • envsupport@licor.com • www.licor.com

Doc. #980-07079