



## Technical Data Sheet

### DeviceLink™ Connect Remote Monitoring Solution Kits

Thermo Fisher Scientific, Asheville, North Carolina

Revision A

Specifications	Catalog Number	
	<b>DLCKTULTS101</b> <b>DLCKTLRFZ101</b> <b>DLCKTCO21101</b> <b>DLCKTCO21102</b>	
<b>Application and Operating Conditions</b>		
Kit Contents	<ul style="list-style-type: none"> <li>• 1x DeviceLink Connect mote (PN: <b>DLCPMTE001</b>)</li> <li>• 1x PoE power supply adapter and PoE cable (PN: <b>DLCPWPWC001</b>)</li> <li>• 1x USB commissioner dongle (PN: <b>DLCSWCOMM001</b>)</li> <li>• 1x Serial/USB cable (PN: <b>Kit dependent</b>)</li> </ul>	<ul style="list-style-type: none"> <li>• 1x Wire cabling/mounting kit (PN: <b>DLCPMONT001</b>)</li> <li>• 1x Quick start installation guide (PN: <b>DLCPINST001</b>)</li> <li>• 1x Year 1 Cloud subscription (PN: <b>DLCSBYEAR001</b>)</li> </ul>
Application	Wireless monitoring solution for laboratory equipment utilizing the Thermo Connect cloud environment for 24/7 remote monitoring and notifications. Direct connection to equipment via serial/USB cable and/or independent sensors	
Application Environment	Non-corrosive, non-flammable, non-explosive, good air ventilation	
Ambient Operating Temperature	Temperature range for exposure of the mote (based on internal component design) Operating temperature range: 0°C to +45°C Operating humidity range: 0% to 95% RH, non-condensing	
Indoor/Outdoor Usage	Indoor use only	
Device Mounting	Device can be mounted to any flat surface. It is recommended to mount to the top of the instrument if possible	
Serial Based Connections	DeviceLink Connect is intended for a 1-to-1 connection with an instrument. Only one serial port based connection per DeviceLink Connect. (Independent sensors can be used in tandem with a serial based connection on a single DeviceLink Connect.)	
<b>Power Ratings and Specifications</b>		
Power Adapter Model	POE16R-1AFG-R	
Power Adapter Rating/Specifications	DC voltage: 56VDC; Load: 0A (min) / 0.275A (max) Regulation Line/Load: 4% Minimum supply power: 15.4W The power supply must conform to the PoE (Power over Ethernet) IEEE802.3af standard	
Power Adapter Wire Length	14 ft (4.27m)	
POE Wall Adapter Dimensions	3.15" x 1.77" x 1.29" (80 mm x 45 mm x 33 mm) (LxWxH)	
POE Wall Adapter Mass	3.35 oz (95 g)	
Backup Battery	2600mAh, 3.6V Rechargeable Li-ion battery providing up to 3hrs of backup power during continuous operation	
<b>Networking Requirements</b>		
Connection Type	Wi-Fi or Ethernet (LAN) connection	
Supported Wi-Fi Protocols	WPA Enterprise (802.1x support); WPA2-PSK	
Wi-Fi Bands and Frequency	Wi-Fi (802.11n/ac) by default; 2.4GHz and 5.0GHz ISM band operation	
Hosting Protocols	DHCP and DNS; NTP local or externally sourced. Device transmits via HTTP(S) periodically per customer configuration.	
Required Ports	Ports 80 (HTTP) or Port 443 (HTTPS) full access to api.traxxekg.com. An explicit HTTP proxy is configurable.	
Supported Web Application Browsers	Internet Explorer™ (10 and later), Google Chrome™ (25 and later), Mozilla Firefox™ (16 and later), Apple Safari™ (7 and later)	
<b>Mote Hardware Design and Connection Ports</b>		
Main Controller	Raspberry Pi 3B+ module	
Indicators and Buttons	LED Indicators (3bi-color red/green): Power/Status, Network, and Sensor Action Button: Manually send sensor data to mote Power Switch: Power mote On/Off	
LAN Ports	LAN (Label: LAN): 10/100 Ethernet (Used for device commissioning only)	
USB Ports	LAN (Label: POE LAN): 10/100 Ethernet with PoE (used for device power and LAN network connection)	
RS485/232 Port	4x USB ports for sensor and cable connection	
Dimensions	1x port for digital inputs (door switches and some sensors)	
Mass	4.82" x 3.78" x 1.50" (122 mm x 96 mm x 38 mm) (L x W x H)	
	10.4 oz (295 g)	
<b>Datalogging, Alarms, and Notifications</b>		
Device Storage	8 GB; Approximately 7 days storage capacity during a network outage <sup>1</sup>	
Remote Access Platform	InstrumentConnect™ mobile and web platforms	
Supported Web Browsers	Internet Explorer™ (10 and later), Google Chrome™ (25 and later), Mozilla Firefox™ (16 and later), Apple Safari™ (7 and later)	
Notifications	Sent via email, In-application notification, and mobile push notifications	
Reports	Reports in either .csv or .pdf available for download via InstrumentConnect web-based application	
<b>Standards and Regulations</b>		
FCC / IC	Contains FCC ID: 2ABCB-RPI3BP; Contains IC: 20953-RPI3P	
EMC	Compliance per Declaration of Conformity 47 CFR § 2.1077 FCC 47 CFR Part 15B, ISED ICES-003	
Safety	UL61010-1:2012, CSA C22.2 No. 61010-1: 2012 from TUV SUD	
Power Adapter Agency Listing	Fully compliant with IEEE802.3af	
Country of Origin	Assembled in USA	

<sup>1</sup> Internal datalogging capacity during a network outage depends on the number of connected sensors and the frequency of data transmission from those sensors.

For a list of serial-based connection compatible equipment visit [thermofisher.com/DLCEquipment](http://thermofisher.com/DLCEquipment)

© 2022 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.