

InfinityLab ERI Remote and Leak Module for connections of up to 4 external leak sensors and one APG remote cable

Installation

This Technical Note describes the installation of the 5067-6149 *External Leak Sensor Assembly* and the 5188-8053 *ERI Leak and Remote Assembly*, to an Agilent Infinity II system. The External Leak Sensor Assembly allows the Agilent Infinity II system to be shut down, and flow to stopped, if a leak has been detected away from the Agilent Infinity II system, such as near large columns or with 3rd party modules.

The ERI Leak and Remote Assembly allows the APG-Remote function of the ERI (Enhanced Remote Interface) Connector of a module to be extended while increasing the number of Leak Sensor ports.

External Leak Sensor Assembly



Figure 1 External Leak Sensor Assembly

Item	p/n	Description
1	5067-6149	Leak Sensor Assembly, 3 m
2		ERI Remote port, male
3		ERI Remote port, female
4		Leak Sensor port (4x)

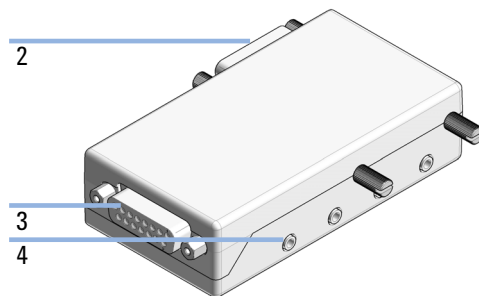
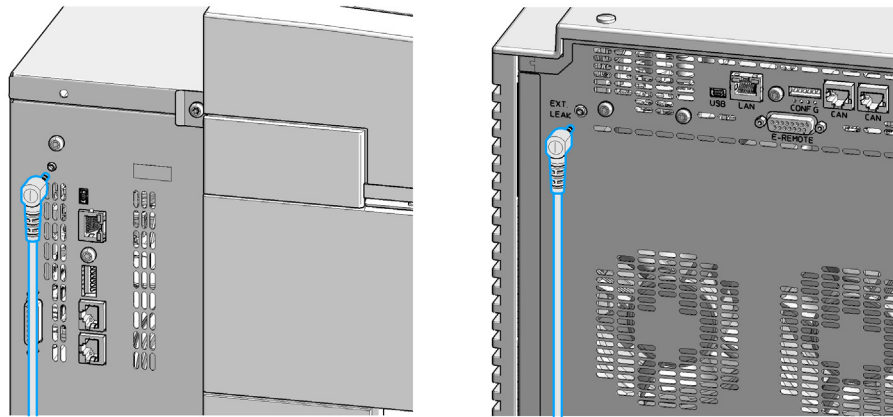


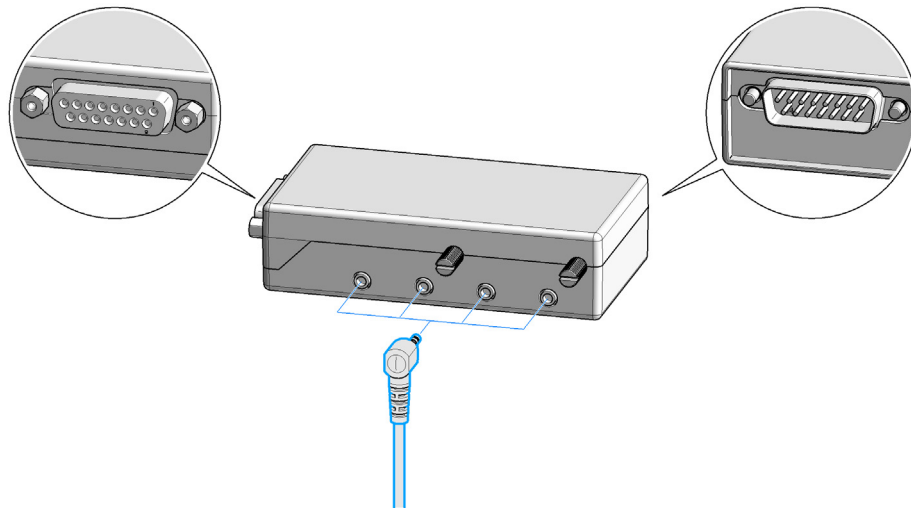
Figure 2 ERI Leak and Remote Assembly

Install the External Leak Sensor Assembly and ERI Leak and Remote Assembly

- 1 Identify the External Leak Sensor Connector on the side of G7158 or G7159, or on the back of the G7161 module, as to where to install the External Leak Sensor assembly, and insert the Connector.



- 2 Alternatively, you can connect a Leak Sensor to the ERI Remote / Leak Module (PN 5188-8053), which allows you to connect up to 4 External Leak Sensors. In addition the ERI Remote / Leak Module offers an extension of the APG Remote function via a 15 Pin Connector.



NOTE

After connecting the External Leak Sensor, wait about 1 min to allow the Leak Sensor to be recognized.

