

Identify the time to change the filter

The Gas Clean sensor LED changes from green to yellow when it is time to change the filter.

	Oxygen filter	Moisture filter/ Process moisture filter	Carrier gas filter	Sulfur filter
	Removes oxygen as well as traces of sulfur and chlorine compounds from carrier gas	Removes water, oil, and other foreign material from carrier gas	Single combination filter. Removes water, oxygen, and organic compounds	Removes water and sulfur-specific compounds such as H ₂ S, COS, and SO ₂ from gas streams
Indicator color change	From green to gray	From green to brown	Oxygen: green to gray Moisture: green to pale brown	From green to pale brown

© Agilent Technologies, Inc. 2019
Printed in the USA, January 2, 2019

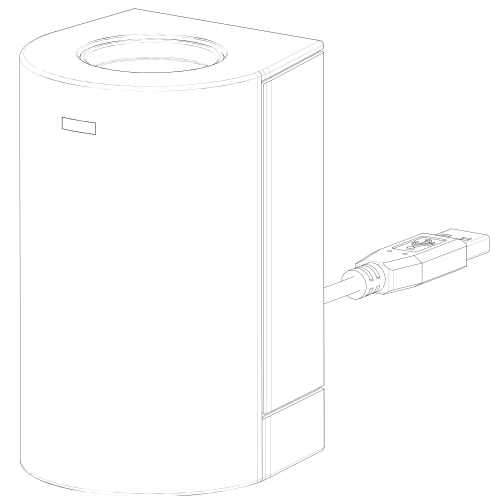


A Step-by-Step Guide to Installing the Gas Clean Sensor

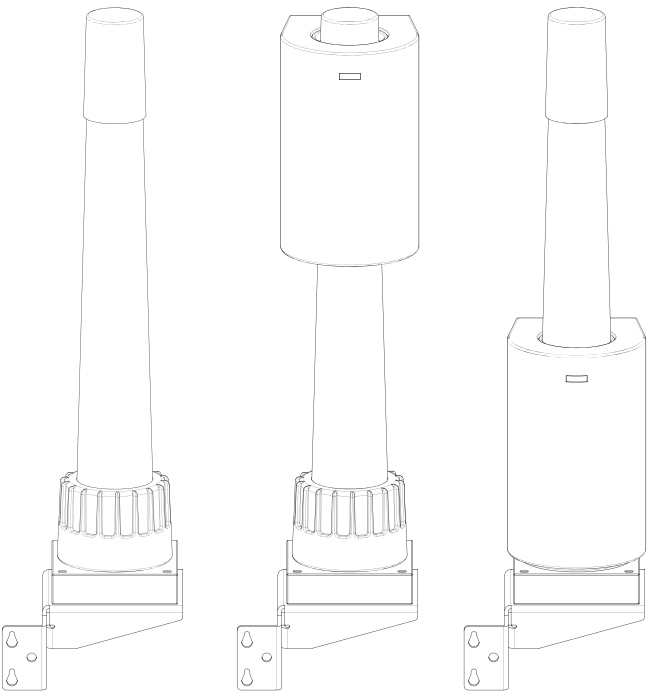


Installing the Gas Clean filter and bracket onto the GC

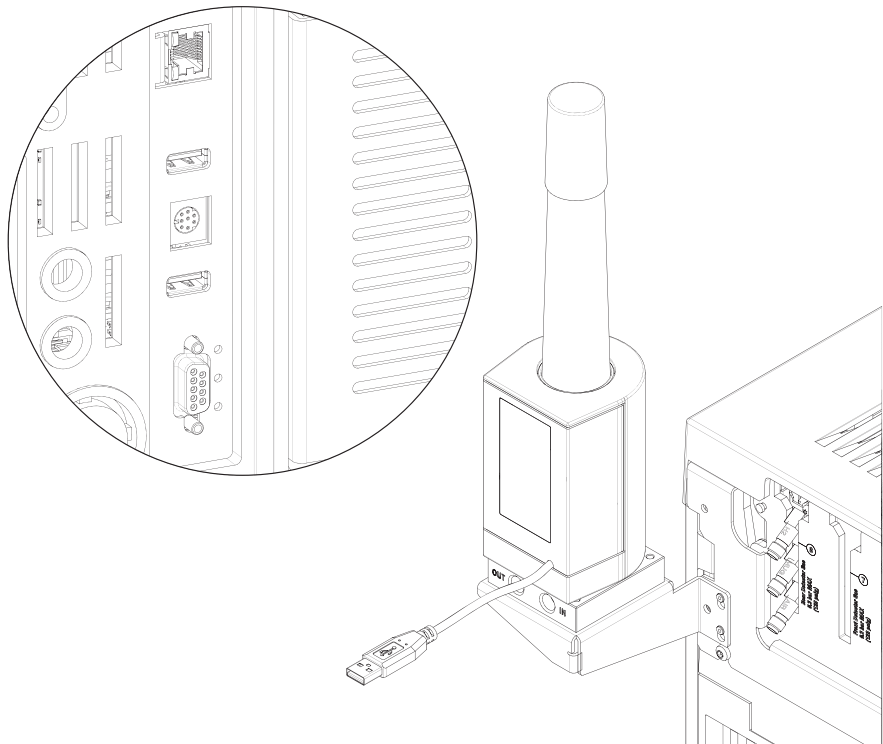
Step 1: Remove the Gas Clean sensor from its packaging.



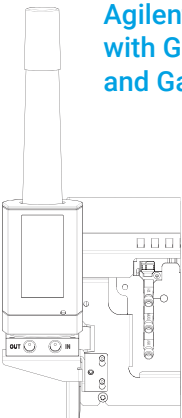
Step 2: Slide the Gas Clean sensor onto the Gas Clean filter.



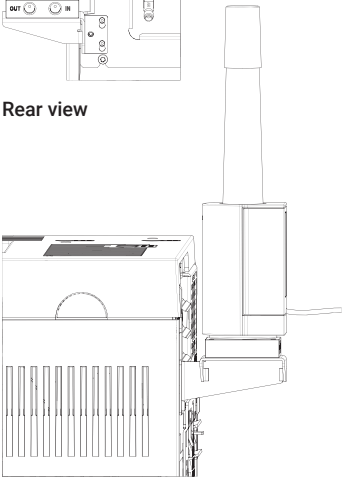
Step 3: Plug the USB cable from the Gas Clean sensor into the USB port.



Agilent 8890 GC
with Gas Clean filter
and Gas Clean sensor



Rear view



Front view