

To Configure GC-HS Communications

The headspace sampler can be configured to communicate with the connected GC. This feature requires a HS with firmware A.01.06 (or greater) and a compatible Agilent GC (for example, a 7890B GC with firmware B.02.02 or greater). With communications configured, the HS knows GC method timings and programs, synchronizes with the GC clock, and follows the GC instrument schedule.

To configure GC-HS communications:

1. At the headspace sampler keyboard, press [**Options**], then go to **Communications > GC**.
2. Enter the GC's IP address, then press [**Enter**].
3. At the GC keyboard, press [**Front Injector**] if the headspace sampler is connected to the GC front inlet, or press [**Back Injector**] if it is connected to the GC back inlet.
4. Scroll to **No Headspace**, then press [**Mode/Type**]. A list of headspace sampler models appears. Scroll to the correct one, then press [**Enter**].
5. Press [**Options**], then select **Communications**.
6. Scroll to the IP address entry for the headspace sampler, either **Fr. Headspace** or **Bk. Headspace**. Enter the headspace sampler's IP address.
7. Press [**Front Injector**] or [**Back Injector**] as applicable.
8. A new **Communication** line appears. Scroll to it and press [**On/Yes**]. The line changes to read **Connected time**. Scroll down to display a read-only summary of the headspace sampler configuration which the GC obtained from the sampler.

If communications fails, check the IP addresses. Verify that the GC's IP address is correct in the headspace sampler, and that the headspace sampler's IP address is correct in the GC. Also check that both instruments are turned on and connected to the LAN. Verify all LAN cable connections.

To temporarily disable communications:

1. At the GC keyboard, press [**Front Injector**] or [**Back Injector**] as applicable.
2. Scroll to **Connected time** and press [**Off/No**].

This setting disables communications but leaves the configuration information intact.