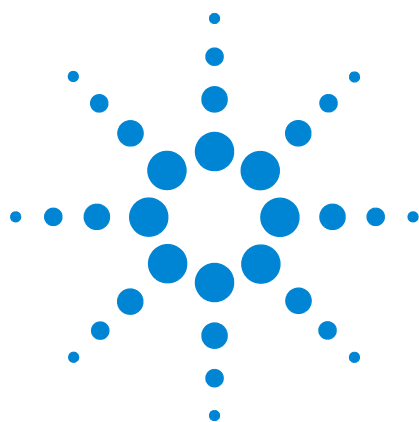




Agilent Multi Mode Inlet Cleaning Guide

Table 1 Parts

Part number	Description	Quantity
G3510-80821	Abrasive brush	1
N/A	Acetone or methanol	25 mL
N/A	100 mL beaker	1

Cleaning Procedure

This procedure explains how to clean the Multi Mode Inlet (MMI). Depending on the inlet mode used, the liner installed, and the cleanliness of the sample, the frequency of cleaning may range from weekly to yearly. When establishing your cleaning frequency, start with a visual inspection of the inlet bottom whenever a liner is changed. A small ring of material collects at the bottom when dirty samples such as food extracts or solid waste extracts are injected. An initial cleaning schedule of every two weeks for dirty samples and every two months for clean samples is appropriate and can be adjusted subsequently.

- 1 Cool the oven and MMI. Turn off the GC for safety.
- 2 Open the MMI turn top assembly and remove the liner. From inside the oven, uninstall the column from the bottom of the MMI.
- 3 Moisten an abrasive brush with solvent and wipe the inside bottom seal of the inlet repeatedly to remove any loose material from the MMI. Rinse the brush with clean solvent to remove any loosened material.
- 4 Place the brush all the way to the bottom of the inlet and, with moderate pressure, twist the brush clockwise and counter-clockwise while it is in contact with the bottom of the MMI (approximately 20 cycles should be sufficient).

NOTE

The abrasive brush can be used multiple times.





- 5** Rinse the brush with clean solvent. Depending on the amount of use and maximum temperature, the bottom of the inlet may still show some oxidation (light brown color). This is normal and will not affect inlet performance. Repeat steps 3–5 (if necessary) to remove any remaining deposits.
- 6** Place a beaker or other container underneath the outlet of the MMI. Using clean solvent, rinse the inlet bore several times but never allow the liquid to fill more than half of the bore. Catch the used solvent at the bottom of the MMI and dispose of it in an appropriate manner.
- 7** Allow the MMI to dry before reinstalling the column and inlet liner.

