



Guard Column

NOTICE: This document contains references to Varian. Please note that Varian, Inc. is now part of Agilent Technologies. For more information, go to www.agilent.com/chem.



This guard column is designed to be used in conjunction with the relevant analytical columns. It should be connected immediately prior to the first analytical column to provide protection against blockage by particulate matter and other damaging material in the eluent or sample. It is recommended that eluents and samples are filtered (0.5 μm) prior to use. The same flow rate, pressure and solvent restrictions detailed in the Users Guide apply to Guard Columns.

PTO for Diagram

SI-02376/A/1.10 Printed in the UK

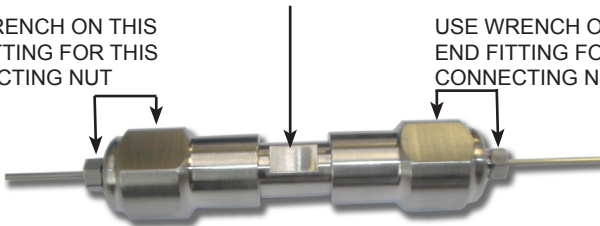




DO NOT USE WRENCH
ON THIS FLAT

USE WRENCH ON THIS
END FITTING FOR THIS
CONNECTING NUT

USE WRENCH ON THIS
END FITTING FOR THIS
CONNECTING NUT



WARNING

To avoid loosening end fittings and causing column leakage, please use wrenches as indicated above when connection the guard column to the system.

PTO for Instructions for Use

