

Agilent MetaCarb H Plus
Guard Column
H+ Form
Guard Column Care Information
Part No. A5216

Precautions:

Mobile phase: Use only aqueous acid mobile phases. AVOID ORGANIC SOLVENTS. Samples may contain small amounts organic solvents. If there are any questions, please do not hesitate to call for more information.

pH range: 1-7

Pressure limit: Do not exceed the recommended back pressure on the analytical column with the guard column in place. Prolonged high pressure and/or flow rates may cause a void in the guard column polymer bed. Replacement of the Agilent A5211 Guard column may become necessary if the guard column back pressure is too high.

Maximum mobile phase flow rate: Do not exceed recommended analytical column flow rate.

Storage solvent:

Guard columns may be stored in mobile phase overnight. Long term storage solvent is in $0.001\ N\ H_2SO_4$. Storage in mobile phase may have corrosive effects on the guard column body and may lead to reduced capacity and/or high column backpressure. DO NOT LET THE GUARD COLUMN DRY OUT. Use guard column end plugs.



Using the guard column:

Mobile phases: Mobile phase of choice is dilute sulphuric at a concentration between 0.0001 and 0.05 N.

Regeneration procedures:

Anion or Metal contamination: Guard column contamination is characterized by high backpressures. Pump the guard column (inverted) with 10 mL of 0.05 N H_2SO_4 . Severe contamination will require pumping in the back flush mode overnight at 0.5 mL/min. Re-equilibrate the guard column with the mobile phase.

Organic contamination: Use a Guard-disc to prevent contamination. To regenerate, pump $0.1\ N\ H_2SO_4$ at $0.5\ mL/min$ overnight.

Bacterial growth: No regeneration procedure is available. To prevent bacterial growth, store the guard column in acidic solution and filter samples and mobile phase through 0.2 µm membrane.

www.agilent.com

Agilent shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance, or use of this material

Information, descriptions, and specifications in this publication are subject to change without notice.

©Agilent Technologies, Inc. 2013

A5216

5991-3017EN

