



**Agilent Technologies**

## **GC Syringe selection Guide.**

This document is believed to be accurate and up-to-date. However, Agilent Technologies, Inc. cannot assume responsibility for the use of this material.  
The information contained herein is intended for use by informed individuals who can and must determine its fitness for their purpose.

