



Agilent 490 Micro GC Natural Gas Analyzer Installation and Configuration

This document describes where to find detailed software installation and configuration information for the Agilent 490 Micro GC Natural Gas Analyzer. See the section(s) below as appropriate for the available software.

Agilent OpenLAB CDS EZChrom Edition or ChemStation Edition

For installation and configuration information, see the *Agilent OpenLAB Chromatography Data System (CDS) Workstation Installation Guide* (CDS_WS-Installation Guide.pdf). Find this document on the OpenLAB CDS DVD in folder **\Disk1\Docs**. Available in several languages.

Agilent EZChrom Elite

For installation and configuration information, see the *EZChrom Elite Installation Guide*.

EZReporter

EZReporter is an optional software package for calculating the calorific value of natural gas samples (heating value calculation). Installation of EZReporter will be done by an Agilent Customer Engineer.

Find detailed information about EZReporter installation and configuration in the following manuals:

- *EZReporter Quick Start Guide* (either the *EZChrom Edition* or *ChemStation Edition*)
- *EZReporter Software* reference manual





To establish the connection between EZReporter and the Agilent data system, two steps are important:

- 1 Calibration.** For creating a calibration method or performing calibration runs, find detailed instructions in either:
 - *EZChrom Elite online help and tutorial*
 - *OpenLAB CDS online help and tutorial*
- 2 Exporting Raw Data.** Configure the Agilent data system (EZChrom Elite or OpenLAB CDS, either edition) to export raw data as described in:
 - *EZReporter Quick Start Guide—EZChrom Edition*
 - *EZReporter Quick Start Guide—ChemStation Edition*

The material contained in this document is provided “as is,” and is subject to being changed, without notice, in future editions. Further, to the maximum extent permitted by applicable law, Agilent disclaims all warranties, either express or implied, with regard to this manual and any information contained herein, including but not limited to the implied warranties of merchantability and fitness for a particular purpose. Agilent shall not be liable for errors or for incidental or consequential damages in connection with the furnishing, use, or performance of this document or of any information contained herein. Should Agilent and the user have a separate written agreement with warranty terms covering the material in this document that conflict with these terms, the warranty terms in the separate agreement shall control.

© Agilent Technologies, Inc. 2012

No part of this manual may be reproduced in any form or by any means (including electronic storage and retrieval or translation into a foreign language) without prior agreement and written consent from Agilent Technologies, Inc. as governed by United States and international copyright laws.

G3582-90003

First edition, May 2012

Printed in The Netherlands

P.O. Box 8033
4330 EA Middelburg
The Netherlands