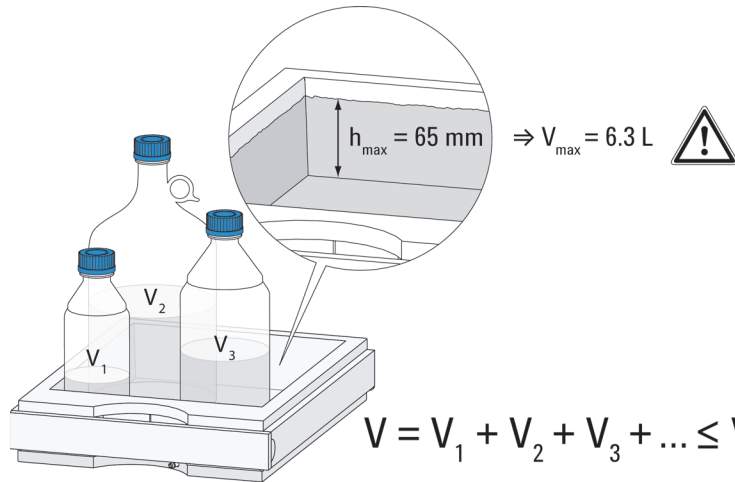
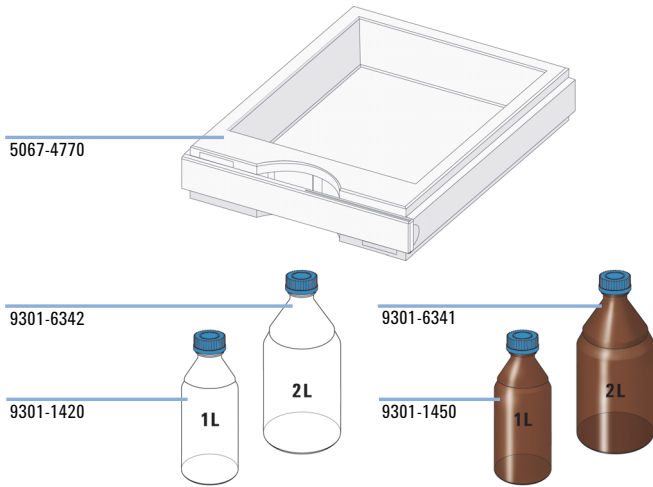
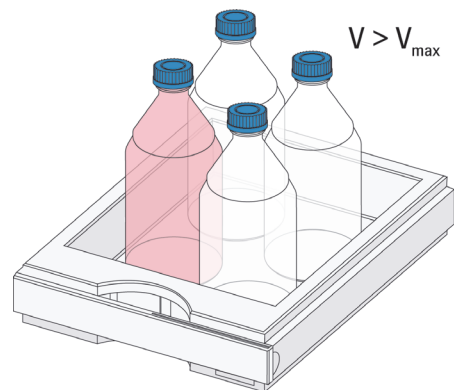
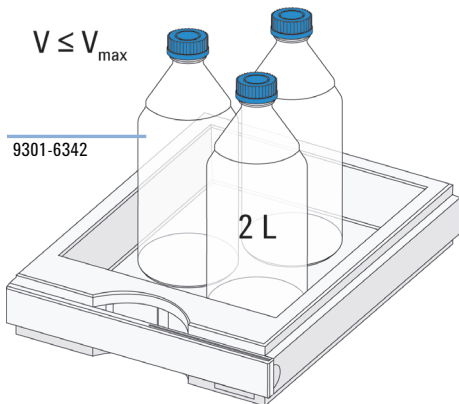
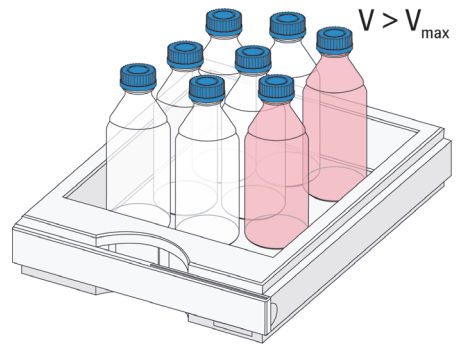
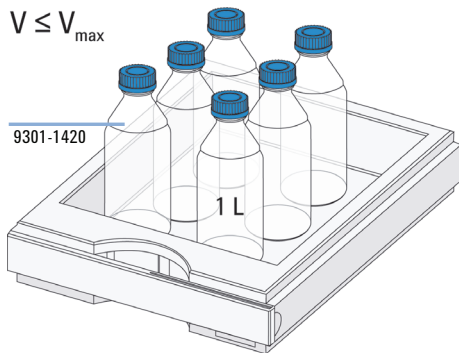
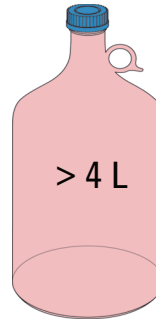
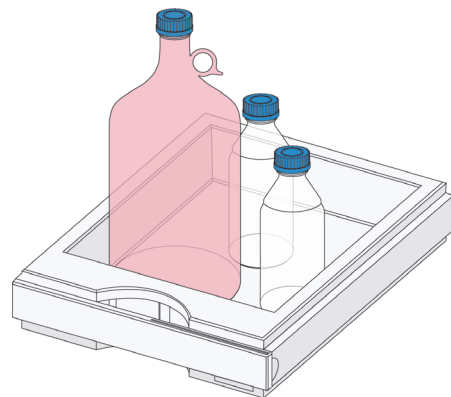
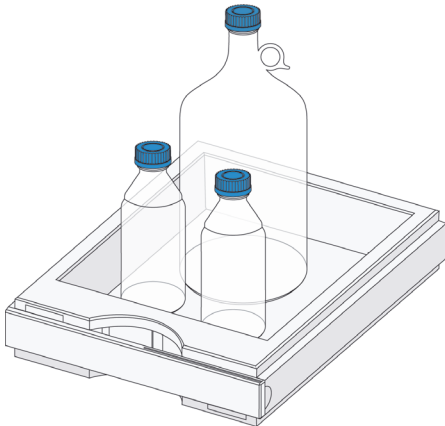
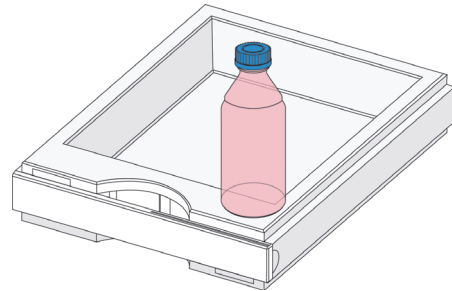
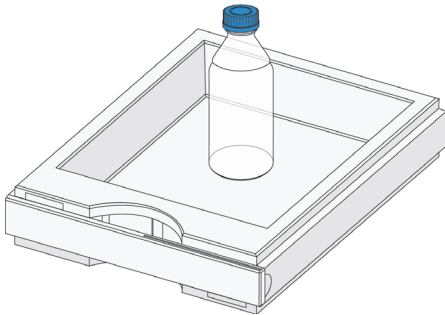
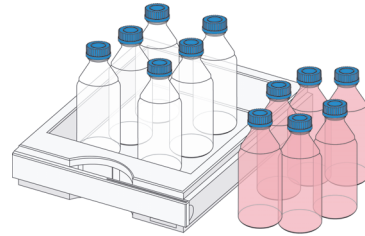
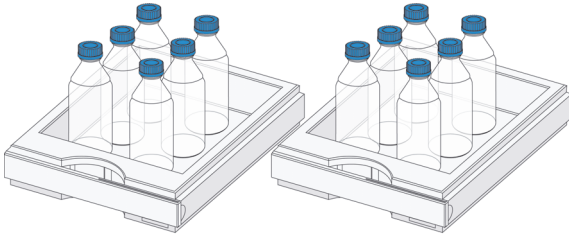


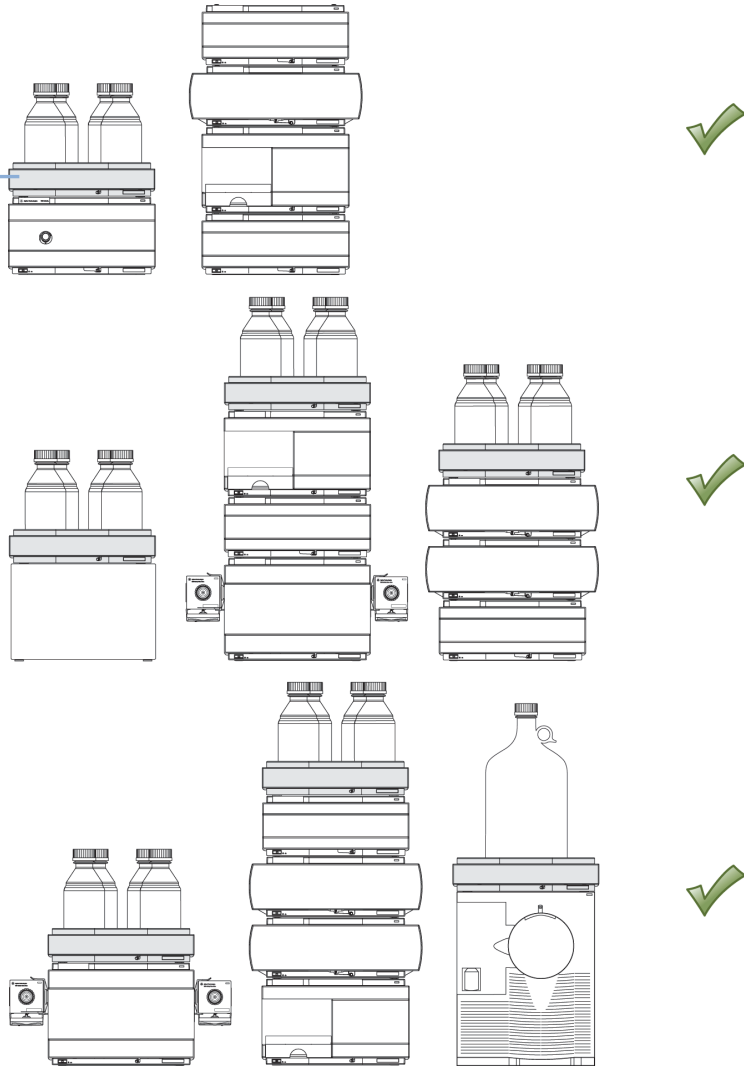
# Usage Guideline for the Solvent Cabinet







5067-4770



01200-90050

Part Number: 01200-90050

Edition 08/2013

Printed in Germany

© Agilent Technologies, Inc. 2013

Agilent Technologies

Hewlett-Packard-Straße 8

76337 Waldbronn, Germany