



Agilent Technologies

UV Cylindrical Cells

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.



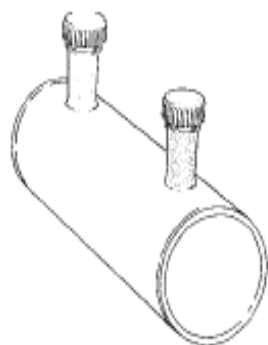
Agilent Technologies

Title: Cylindrical cells

A cylindrical cell is a cell with plane-parallel optical surfaces whose inner volume is cylindrical in shape and has a longitudinal axis parallel to the direction of the radiation beam.

The cells are equipped with two fittings to facilitate filling and emptying of the cells. It is recommended not to remove the cylindrical cell between blank and sample measurement.

Cylindrical Cells with PTFE stopper



Path Length [mm]	External Dimensions L x Diameter [mm]	Internal Diameter [mm]	Volume [μ l]	Part No. Glass	Part No. Quartz
100	102.5 x 22	19	28000	5063-6566	5061-3392

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.