



Agilent Technologies

UV Macro 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: Macro cells

The macro cell, which is defined by DIN 58963 as a rectangular cell with an inner width greater than 5 mm, has emerged as the standard for photometry. The most widely used macro cell is a rectangular cell with outer dimensions of 45 x 12.5 mm (height x width). The length of the cell is dependent on the desired path length. Please refer below for some of the available Macro cells.

Macro Cells with PTFE lid



Path Length [mm]	External Dimensions H x W x D [mm]	Internal Dimensions H x W [mm]	Volume [μl]	Part No. Glass	Part No. Quartz
1	45 x 12.5 x 3.5	44.5 x 9.5	350	5063-6546	5061-3384
2	45 x 12.5 x 4.5	44.5 x 9.5	700	5063-6547	5061-3385
5	45 x 12.5 x 7.5	44.5 x 9.5	1750	5063-6548	5061-3386
10	45 x 12.5 x 12.5	44.5 x 9.5	3500	5063-6549	5061-3387
20	45 x 12.5 x 22.5	44.5 x 9.5	7000	5063-6551	5063-6553
50	45 x 12.5 x 52.5	44.5 x 9.5	17500	5063-6552	5063-6554
Matched pair, set of two:					
10	45 x 12.5 x 12.5	44.5 x 9.5	3500	5063-6550	1000-0544

Macro Cells with PTFE stopper



Path Length [mm]	External Dimensions H x W x D [mm]	Internal Dimensions H x W [mm]	Volume [μl]	Part No. Glass	Part No. Quartz
5	48 x 12.5 x 7.5	42 x 9.5	1750	5063-6555	5063-6557
10	48 x 12.5 x 12.5	42 x 9.5	3500	5063-6556	5062-2477

Note: Spacers are required for cells with outer depth less than 12.5 mm in order to hold them securely in the cell holder.
See page 21.

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.