



Agilent Technologies

Approximate Lamp Lifetime Hours

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

Approximate LAMP Lifetime of UV detectors.

Instrument	Detector	Part number	expected Life time [hours]
HP 1050	79853A VWD	79853-60002	1000
	79853C VWD	79883-60002	1000
	79854A DAD	79883-60002	1000
HP 1090	79880A DAD	79883-60002	1000
	79881A FPD	79880-60002	1000
1100 Series	G1314A VWD	G1314-60100	1000
	G1315A DAD	G1103-60001	2000
	DAD/MWD	2140-0590	1000
	DAD/MWD	5181-1528	2000
	DAD/MWD	5181-1530	2000
G1600A	CE	2140-0585	1000
G1321A	FLD	2140-0600	2000
HP 8452A		08452-60104	1000
8453 Series	Deuterium Lamp	2140-0605	1000
	Tungsten Lamp	G1103-60001	2000

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.