



Agilent Technologies

How to select the Reference Wavelength

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.

How to select Reference Wavelength.

If you know little about the analytes in your sample, use both lamps and store all spectra over the full wavelength range. The detector measures absorbance simultaneously at wavelengths from 190 to 950 nm. Two lamps provide good sensitivity over the whole wavelength range. The deuterium discharge lamp provides the energy for the UV range (190 to 400 nm) and the tungsten lamp emits light from 400 to 950 nm for the visible and short wave near infrared.

Select the Sample Wavelength at or near apex of the peak and select wavelength band with the width of sample bandwidth (BW)

Select the Reference Wavelength in a place in your spectra where your sample is not having any absorbance or zero absorbance and select wavelength band with the width of Reference bandwidth (BW)

Please refer to the Spectra of Anisic acid and use this Spectral information to optimize wavelength settings for your chromatographic signal.

