

# Universal ATR Polarization Accessory for Spectrum Series FT-IR Spectrometers



Attenuated total reflectance (ATR) techniques have been well established in FT-IR spectroscopy for many years. They offer a convenient alternative to thin-film, KBr pellet or diffuse reflectance measurements.

Horizontal ATR accessories using crystal materials such as zinc selenide, thallium bromide/iodide and germanium are used extensively for analysis of liquids, oils, waxes and soft polymeric materials. However, since these crystal materials are relatively soft, they can be easily damaged and are therefore not ideal for the analysis of powders or hard solids.

The introduction of diamond ATR units has meant that the vast majority of sample types, whether liquids, powders, or hard minerals, can be analyzed routinely using a single accessory.

## Key Features

- ▶ Straightforward analysis of most sample types with a minimum of sample preparation
- ▶ Interactive pressure control ensures optimum contact between sample and diamond for high quality, reproducible spectra
- ▶ ATR automatically recognized and checked by instrument
- ▶ Easy to clean diamond top-plate

The PerkinElmer® Universal ATR Accessory offers straightforward analysis of most sample types with a minimum of sample preparation. It features an interactive pressure control so that exactly the right contact is made between the sample and the diamond, resulting in high quality, reproducible spectra.

1. The Spectrum™ 100 or Spectrum 400 FT-IR and FT-NIR systems will automatically recognize the accessory, and with user-defined analysis parameters the accessory performance is checked automatically prior to analysis.
2. The diamond top-plate is easy to clean using a dry or solvent-containing tissue.

The Universal ATR Polarization Accessory combines the above sampling benefits with the ability to acquire polarized IR spectra.

A variable-angle polarizer is slotted into the accessory and the user simply sets the required angle. For non-polarized measurements the polarizer is removed.

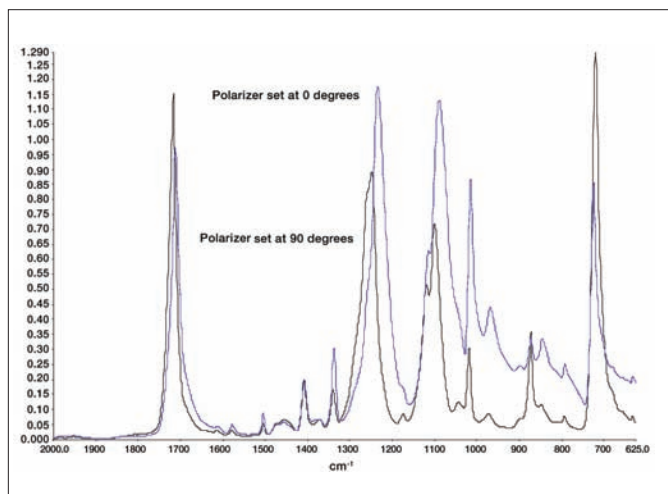


Figure 1. Differences in spectra shown when parallel and perpendicularly polarized spectra from the wall of a PET bottle are measured.

The benefit of polarization measurements is that they give additional information on the molecular structure of a sample, particularly the crystal or molecular orientation. For example, the dichroic ratio is often measured for specific absorbance bands in the spectrum and is defined as the ratio of the absorbance measured using parallel-polarized radiation, divided by the absorbance measured using perpendicular-polarized radiation.

Application areas where these polarization measurements are particularly useful are crystal structure analysis, polymer and fiber orientation studies and structural studies of biomolecules such as proteins.

An example of the need for polarized measurements is in the manufacture of PET (polyethylene terephthalate) bottles. These bottles are blow molded from a pre-form and the strength of the bottle is dependent on the long chain molecules of the PET polymer being preferentially oriented around the circumference rather than along the axis. The polarized IR spectra from the wall of a bottle are shown in Figure 1, demonstrating the difference in the spectra when parallel and perpendicular polarized spectra are measured.

## Ordering information

### Universal ATR Polarization Accessories:

Single reflection accessory	L1250064
Three reflection accessory	L1250065
Nine reflection accessory	L1250066

### Standard Universal ATR Accessories:

Single reflection Diamond/ZnSe accessory	L1250050
Three reflection Diamond/ZnSe accessory	L1250051
Nine reflection Diamond/ZnSe accessory	L1250052
Single reflection Diamond/KRS-5 accessory	L1250053
Single reflection Germanium accessory	L1250054

PerkinElmer, Inc.  
940 Winter Street  
Waltham, MA 02451 USA  
Phone: (800) 762-4000 or  
(+1) 203-925-4602  
[www.perkinelmer.com](http://www.perkinelmer.com)



For a complete listing of our global offices, visit [www.perkinelmer.com/lasoffices](http://www.perkinelmer.com/lasoffices)

©2008 PerkinElmer, Inc. All rights reserved. The PerkinElmer logo and design are registered trademarks of PerkinElmer, Inc. Spectrum is a trademark and PerkinElmer is a registered trademark of PerkinElmer, Inc. or its subsidiaries, in the United States and other countries. All other trademarks not owned by PerkinElmer, Inc. or its subsidiaries that are depicted herein are the property of their respective owners. PerkinElmer reserves the right to change this document at any time without notice and disclaims liability for editorial, pictorial or typographical errors.