

**Model 300 Manual**  
**Continuously Variable ATR**  
**0012-01X**  
Version 2.4



P/N 700-0007

**SPECTRA**TECH

A subsidiary of Thermo Optek Corporation,  
a Thermo Electron Company

website: [www.spectra-tech.com](http://www.spectra-tech.com)  
e-mail: [info@spectra-tech.com](mailto:info@spectra-tech.com)

**In USA**  
**Spectra-Tech, Inc.**  
2 Research Drive  
P.O. Box 869  
Shelton, CT. 06484-0869  
Phone: 203-926-8998  
800-843-3847  
FAX: 203-926-8909

**In Japan**  
**ST Japan**  
1-14-10 Kakigara-cho  
Nihonbashi, Chuo-ku  
Tokyo, 103 Japan  
Tel: (033) (666) 2561  
FAX: (033) (666) 2658

# Table of Contents

## General Information

Introduction .....	2
The Manual .....	2
Theory .....	3-4

## Getting Started

Product Description .....	4
Crystal Holder Assembly .....	5-6
Installation & Alignment .....	7-9

## Using the Model 300

Performing an Experiment .....	10-11
--------------------------------	-------

## Appendix A

Changing to a 25mm Crystal .....	12-14
----------------------------------	-------

## Appendix B

Indices of Refraction .....	15
-----------------------------	----

## Appendix C

Properties of Crystal Materials.....	16
--------------------------------------	----

# General Information

---

## Introduction to ATR Spectroscopy

Attenuated Total Internal Reflectance (ATR) Spectroscopy is a versatile and powerful technique for infrared sampling. Since materials are normally analyzed by ATR with either minimal or no sample preparation, ATR is a rapid technique for obtaining the infrared spectrum of a material. Materials which are either too thick or too strongly absorbing to be analyzed by transmission spectroscopy can be routinely analyzed using ATR spectroscopy. ATR is also useful when only the surface of the material is of interest.

---

## The Manual

This manual is designed as a tutorial to guide you through the installation of the Model 300 Continuously Variable ATR and through a typical ATR spectroscopy analysis. If you have any questions, please contact a Spectra-Tech Technical Representative.

---

## Theory

Internal reflection spectroscopy is a common infrared technique in which the infrared radiation is passed through an infrared transmitting crystal of high refractive index, allowing the radiation to reflect in the crystal one or more times. In this way, an evanescent wave penetrates into the sample in contact with the crystal, producing a spectrum of the sample. In the most common application of this technique, ATR (Attenuated Total Reflectance), the sampling surface is held in a vertical orientation, with sample material placed on one or both sides of the ATR crystal.

# General Information

## Title Here

The following equations define the effective angle of incidence for the Model 300. These calculations can be used to determine the best crystal materials for specific applications.

**See Appendix B for a chart of pre-calculated angle of indices.**

$$\text{effective angle of incidence } (\theta) = \theta_{\text{IRA}} - \sin^{-1} \left[ \frac{\sin(\theta_{\text{IRA}} - \theta_{\text{IRE}})}{n_1} \right]$$

$$\# \text{ of reflections} = l/t \cot \theta$$

$$\text{depth of penetration } (d_p) = \frac{\lambda}{2\pi n_1 [\sin^2 \theta - (n_2/n_1)^2]^{1/2}}$$

$$\text{effective pathlength} = N \times d_p$$

$\theta_{\text{IRA}}$  = angle designated on scale

$\theta_{\text{IRE}}$  = crystal face angle

$\lambda$  = wavelength (mm)

$n_1$  = refractive index of crystal

$n_2$  = refractive index of sample

$\theta$  = crystal face angle (degrees)

$l$  = length of crystal

$t$  = thickness of crystal



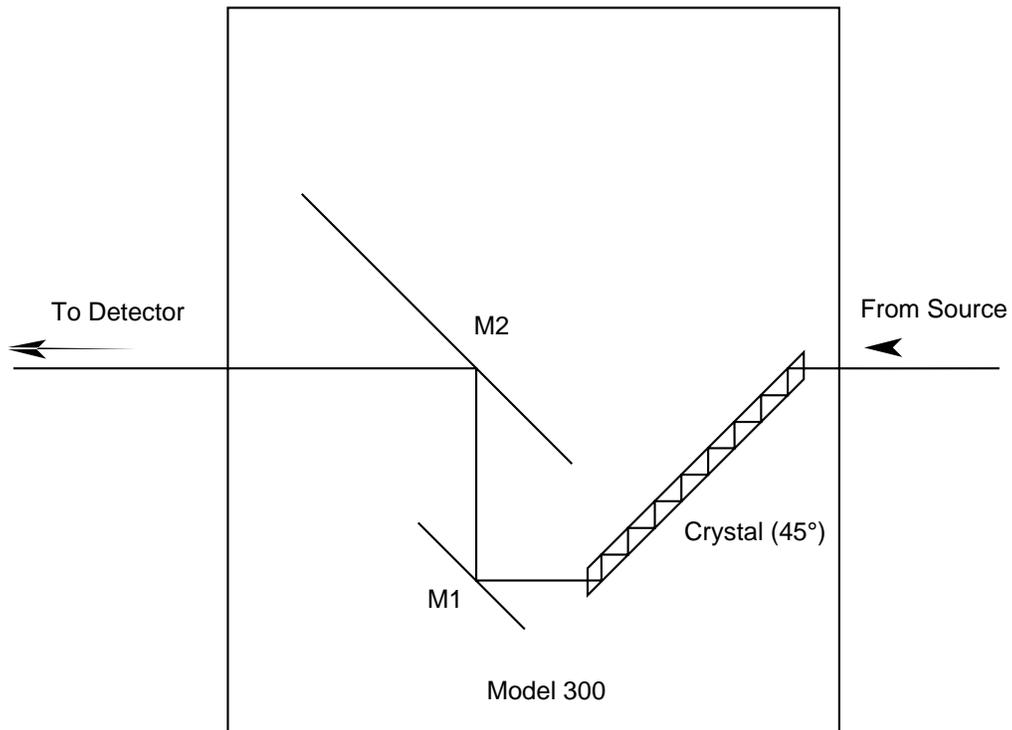
In addition, users will need to verify that the sampling crystal is not soluble or otherwise damaged by the sample. A summary of the physical properties of different crystal materials is listed in Appendix C. If additional information is required contact Spectra-Tech.

For a complete description of the theory of Attenuated Total Reflectance please refer to Spectra-Tech's FT-IR Technical Note #1: *Introduction to Attenuated Total Internal Reflectance (ATR) Spectroscopy*.

# Getting Started

## Product Description

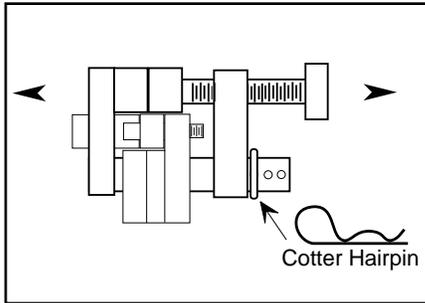
The Model 300 - Continuously Variable ATR is a high performance, continuously variable angle ATR accessory. The angle is continuously variable from  $30^\circ$  to  $60^\circ$ . Unique mirror placement and pivoting make changing angles simple and repeatable while the focus is maintained at all angular positions.



*The diagram above shows the ray path of the IR beam. The energy beam from the source bounces through the crystal, reflects off the surface of mirror #1 onto mirror #2 and out to the detector.*

# Getting Started

## Crystal Holder Assembly

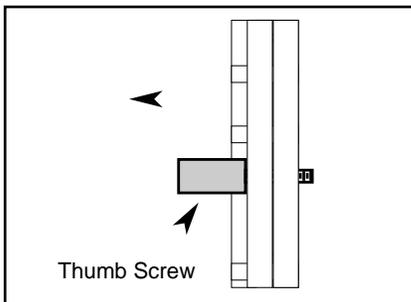


### **Prepare holder for crystal**

Place the sample holder on a flat surface.

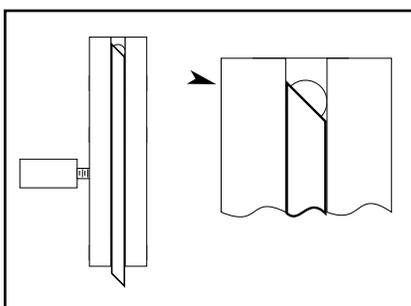
Remove the Cotter Hairpin.

Remove the C-clamp pressure plate assembly and set it aside.



### **Prepare holder for crystal**

Loosen the thumb screw on the crystal holder to open the crystal retention plates.



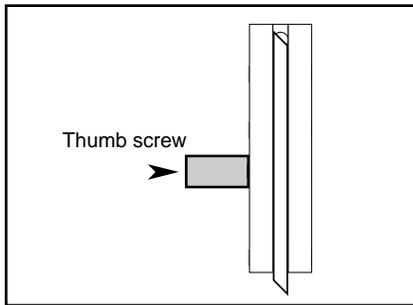
### **Install crystal**

Place the crystal in the holder (in between the retention plates) so that the angled end of the crystal bisects the hole in the base of the holder.

**Note:** Use finger cots to prevent transferring skin oil to the crystal.

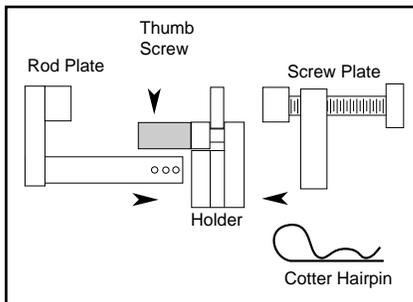
# Getting Started

## Crystal Holder Assembly



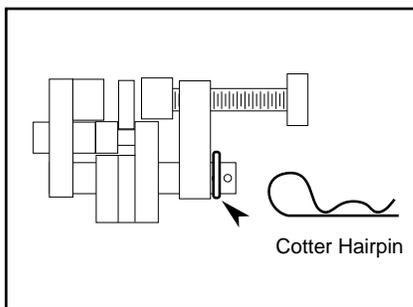
### **Tighten thumb screw**

Finger tighten the thumb screw to hold the crystal in place.



### **Replace C-clamps**

Replace the C-clamp pressure plates- (Place rod plate so that rod fits through hole in holder and thumb screw on holder fits through hole in rod plate. Place screw plate so that rod fits through the hole in the screw plate.)

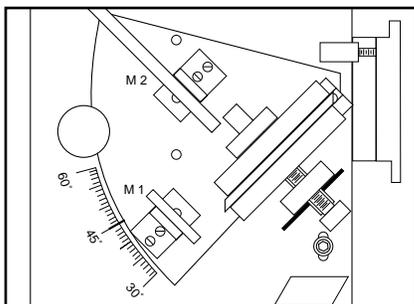


### **Replace locking pin**

Place the Cotter Hair Pin through the hole in the rod.

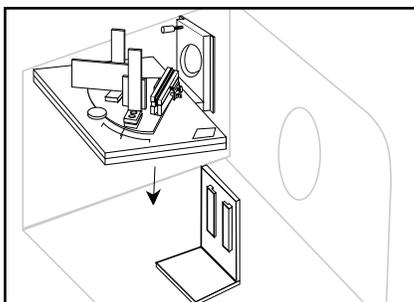
# Getting Started

## Installation & Alignment



### **Place crystal holder in accessory**

Place the crystal holder in the Model 300 so that the hole on the bottom of the holder is sitting on the pin closest to the slide mount and the slot in the bottom of the holder is sitting on the pin closest to mirror #1.



### **Place accessory in instrument**

Place the Model 300 in the sample compartment so that it is sitting securely on the slide mount. Tighten the locking screw (located on the slide mount holder).

### **Set instrument for alignment**

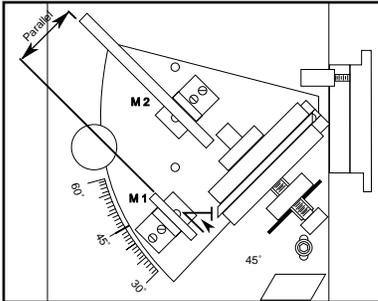
Set your instrument to the energy monitoring mode (e.g. energy, align. tune)

**Note:** If the energy is not greater than 25%\* of the open beam energy, follow the procedure outlined below.

\*With a Ge or KRS-5 crystal, the energy throughput will be lower.

# Getting Started

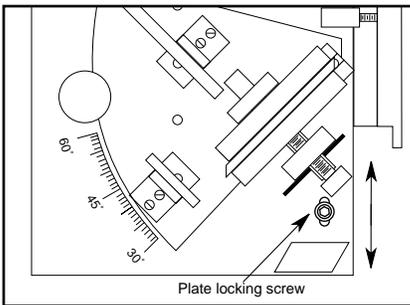
## Installation & Alignment



### **Prepare the accessory for alignment**

Place the angle indicator to 45°.

Adjust the two mirrors so that they are parallel to each other and perpendicular to the crystal.

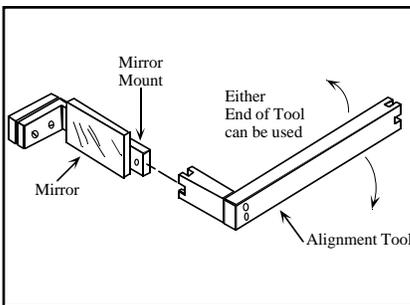


### **Maximize energy**

Using an allen key, loosen the socket head cap screw (SHCS) located above the Spectra-Tech logo.

Slide the accessory's top plate forward and back maximizing the energy.

Tighten the screw.



### **Maximize energy**

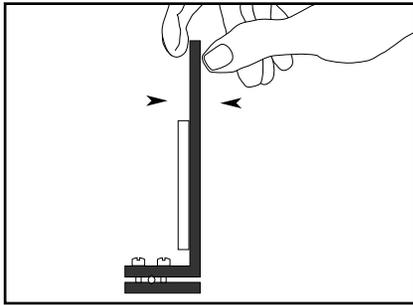
Using the alignment tool, rotate mirror #1 maximizing the energy level.

Using the alignment tool, rotate mirror #2 maximizing the energy level.

Repeat mirror #1 alignment, then mirror #2 alignment.

# Using the Model 300

## Installation & Alignment



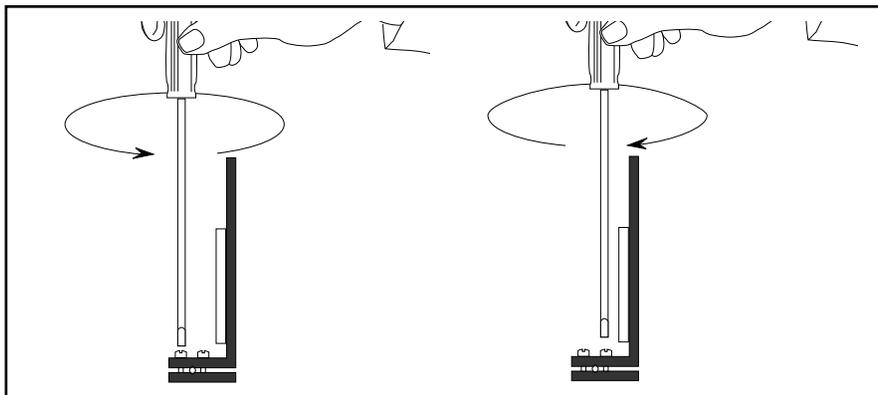
### Adjust the mirror tilt

If the energy level is still below 25%, adjust the tilt on the transfer mirror assemblies as follows: While monitoring either the throughput energy number or the height of the centerburst of the interferogram, gently press down on the front side of the input mirror bracket. Then, repeat this step while pulling gently on the rear side of the mirror bracket.

- If energy increases by applying pressure to the front of the mirror bracket, loosen the front screw and tighten the back screw (see diagram below).
- If energy increases by applying pressure to the back of the mirror bracket, loosen the back screw and tighten the front screw.

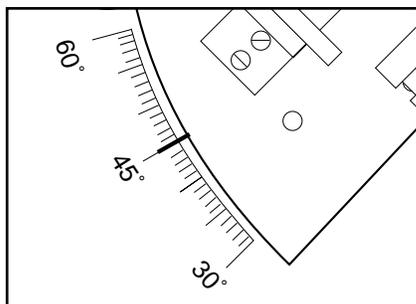
Repeat this step for both transfer mirror assemblies. ***Do not attempt to make this adjustment on just one mirror.***

When both mirror assemblies have been adjusted, check to see that the throughput is greater than 25% of the open beam energy reading.



# Using the Model 300

## Performing an Experiment

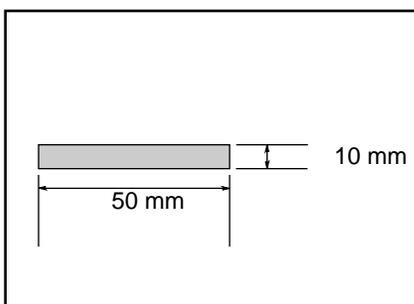


### **Set angle\* & Acquire a background spectrum**

Set the angle indicator to the desired angle setting (45° is the most commonly used setting - see page 15 for calculated angles of indices).

Acquire a background single-beam spectrum with the empty holder in place.

**NOTE:** It is very important to make sure that nothing is touching the crystal faces while the background scan is taken.

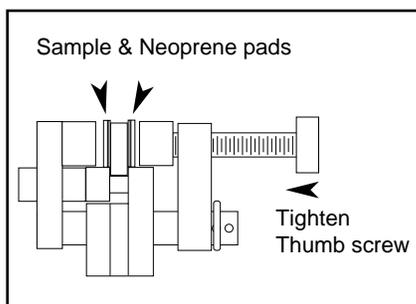


### **Prepare Sample**

Remove the crystal holder from the accessory.

Cut the sample into 2 strips approximately 50mm x 10mm.

**CAUTION:** Be sure that the sample does not interfere with the crystal faces (angled ends).



### **Load Sample**

Place the sample on both sides of the crystal (sandwich the crystal).

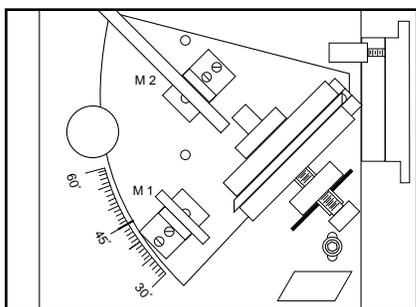
Place the neoprene pads between the sample and the pressure plates.

Tighten the thumb screws to hold the sample in place.

\*For an in depth discussion of angles of incidence and depth of penetration, see Spectra-Tech's Technical Note #1.

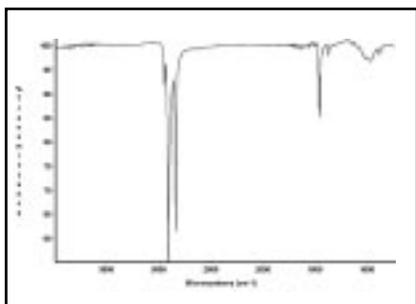
# Using the Model 300

## Performing an Experiment



### **Replace crystal holder**

Replace the crystal holder in the Model 300 so that the hole on the bottom of the holder is sitting on the pin closest to the slide mount and the slot in the bottom of the holder is sitting on the pin closest to mirror #1.

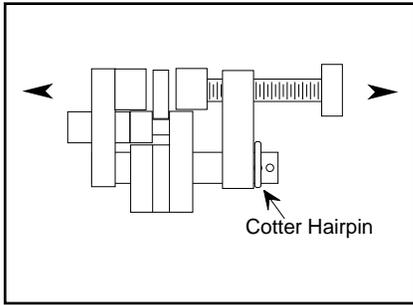


### **Acquire a sample spectrum**

Acquire a sample single-beam spectrum.  
Ratio it against the previously acquired background single-beam spectrum.

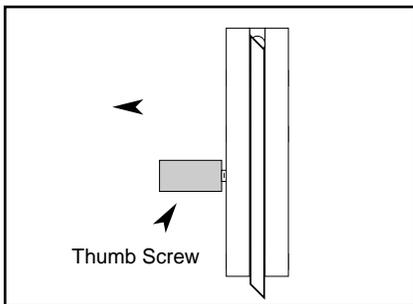
# Appendix A

## Changing to a 25mm crystal



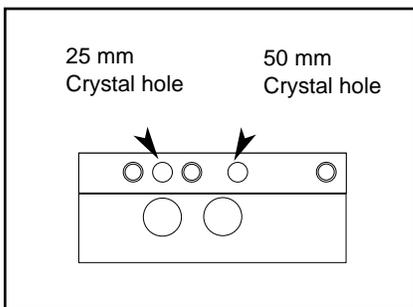
### Prepare holder for crystal

Remove the sample holder from the accessory.  
Place the sample holder on a flat surface.  
Remove the Cotter Hair Pin.  
Remove the C-clamp pressure plate assembly and set it aside.



### Prepare holder for crystal

Loosen the thumb screw on the crystal holder to open the crystal retention plates.  
Remove the 50mm crystal.

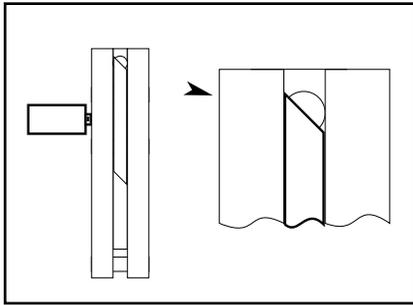


### Prepare holder for crystal

Remove the thumb screw on the crystal holder and shift it to the 25mm location..

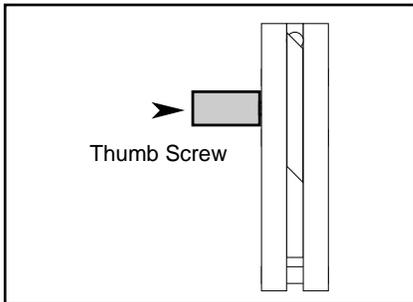
# Using the Model 300

## Changing to a 25mm crystal



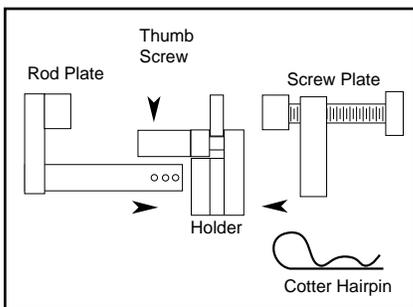
### Install crystal

Place the 25mm crystal in the holder (in between the retention plates) so that the angled end of the crystal bisects the hole in the base of the holder.



### Tighten thumb screw

Tighten the thumb screw to hold the crystal in place.

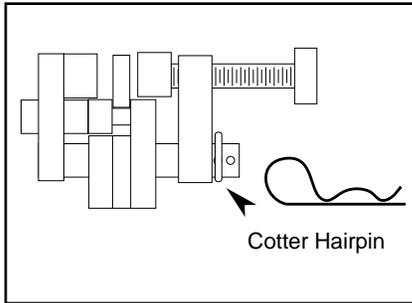


### Replace C-clamps

Replace the C-clamp pressure plates- (Place rod plate so that rod fits through hole in holder and thumb screw on holder fits through hole in rod plate. Place screw plate so that rod fits through the hole in the screw plate.)

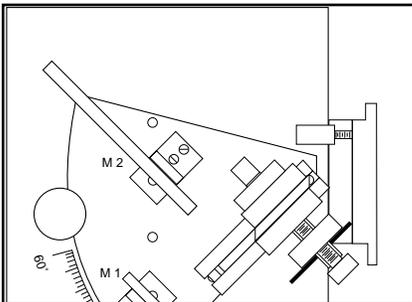
# Using the Model 300

## Changing to a 25mm Crystal



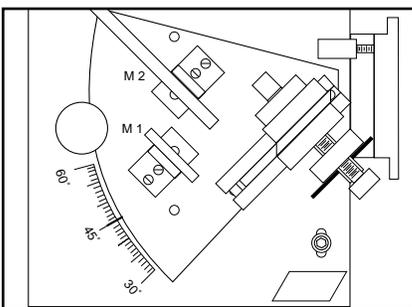
### **Replace locking pin**

Place the Cotter Hair Pin through the hole in the rod.



### **Place crystal holder in accessory**

Place the crystal holder in the Model 300 so that the hole on the bottom of the holder is sitting on the pin closest to the slide mount and the slot in the bottom of the holder is sitting on the pin closest to mirror #1.



### **Change the mirror positions**

Loosen the screw on the base of mirror #1, move mirror #1 to its alternate pin position and tighten the screw.

Follow alignment procedure described on pages 8 and 9.

# Appendix B

## Angles of Incidence allowing for the refractive index of the ATR crystal

$\theta_{IRA}$  = Angle designated on scale

$\theta_{IRE}$  = Calculated Angle of Incidence

**KRS-5** ( $n= 2.38$ )

	<b>30°</b>	<b>45°</b>	<b>60°</b>	$\theta_{IRE}$
$\theta_{IRA}$ <b>30°</b>	30°	38.7°	47.9°	
<b>45°</b>	36.3°	45°	53.7°	
<b>60°</b>	42.1°	51.3°	60°	

**ZnSe** ( $n= 2.40$ )

	<b>30°</b>	<b>45°</b>	<b>60°</b>	$\theta_{IRE}$
$\theta_{IRA}$ <b>30°</b>	30°	38.8°	48°	
<b>45°</b>	36.2°	45°	53.8°	
<b>60°</b>	42°	51.2°	60°	

**Ge** ( $n= 4.01$ )

	<b>30°</b>	<b>45°</b>	<b>60°</b>	$\theta_{IRE}$
$\theta_{IRA}$ <b>30°</b>	30°	41.3°	52.7°	
<b>45°</b>	33.7°	45°	56.3°	
<b>60°</b>	37.3°	48.7°	60°	

**Si** ( $n= 3.40$ )

	<b>30°</b>	<b>45°</b>	<b>60°</b>	$\theta_{IRE}$
$\theta_{IRA}$ <b>30°</b>	30°	40.6°	51.5°	
<b>45°</b>	34.4°	45°	55.6°	
<b>60°</b>	38.5°	49.4°	60°	

# Appendix C

## Properties of Crystal Materials

	ZnSe	Si	Ge	KRS-5
<b>Transmission Range</b> ( $\text{cm}^{-1}$ )	20,000-650	8300-660	5500-600	20,000-250
<b>Refractive index</b> @ $1000 \text{ cm}^{-1}$	2.4	3.4	4.0	2.37
<b>Density</b> ( $\text{g/cm}^3$ )	5.27	2.33	5.32	
<b>Hardness</b> (Knoop #)	150	1150	1150	40.2
<b>Cleaning Agents</b>	acetone, $\text{H}_2\text{O}$	acetone, $\text{H}_2\text{O}$	toluene, $\text{H}_2\text{O}$	MEK
<b>Solvents which attack material</b>	acids, strong alkalies	$\text{HF}+\text{HNO}_3$	hot $\text{H}_2\text{SO}_4$ , aq. regia	complex.agents
<b>Remarks</b>	<i>hard, easily cracked</i>	<i>withstands thermal or mechanical shock</i>	<i>hard and brittle, reflection losses</i>	<i>deforms under pressure, toxic</i>