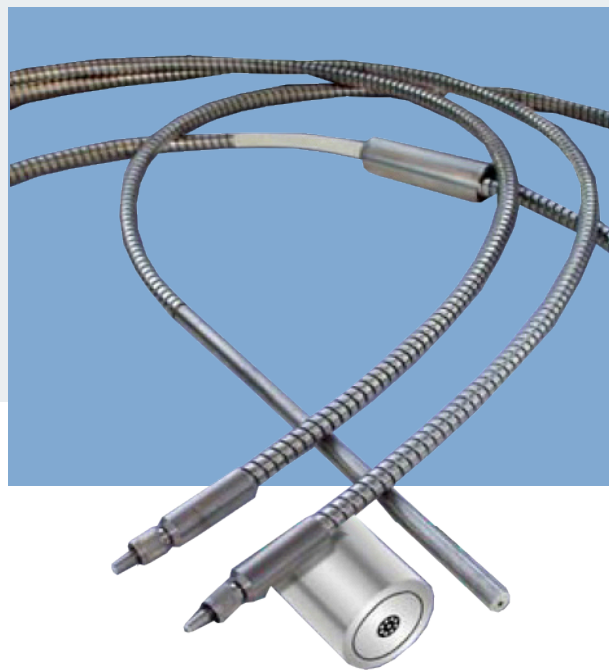


Six-Around-One Reflectance Probe

Spectroscopy



Features

- Economical
- High optical throughput
- Rugged stainless-steel construction
- Small ¼ in. diameter
- SMA fiber connectors
- Compatible with all PerkinElmer spectrometers

The sample is illuminated by six fibers surrounding a single pickup fiber. These probes offer high optical throughput, are rugged, and are easy to use. Standard fiber size is 400 micron core.

Specifications

Part Number: Probe-7-400-SMA

Optical

Numerical aperture 0.22 NA
 Optical fiber type UV or NIR 400/440 silica on silica
 Fiber packaging PVC coated armored duplex
 Fiber length Three-meter standard
 Fiber termination SMA 905

Physical

Body and wand materials 316 SS
 Temperature range -20 °C to 100 °C
 Maximum pressure 250 psi/17 bar

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



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2019, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.