

Spectroscopy

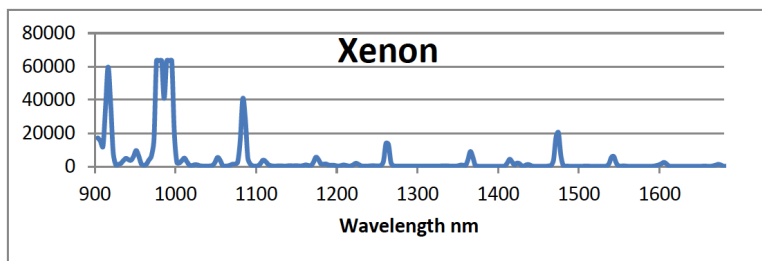
Xenon Calibration Lamp



A calibration source with a diffuse aperture that provides even launch conditions into the fiber-optic cable, resulting in a stable spectrum and sharper lines. This xenon source is ideal for calibrating spectrometers, specifically those operating in the NIR region. A long-life 5000-hour bulb ensures low maintenance. The xenon source can be externally triggered, which is a necessary feature for OEM system integrators.

Xenon Emission Lines (nm)

916.265	1262.339	1672.815
979.970	1365.706	1732.577
992.319	1473.281	2026.224
1083.837	1541.839	



- Diffuse aperture
- Fast, accurate calibrations
- 905 SMA connector
- Trigger
- Regulated power supply
- Compact and rugged

Specifications

Part Number	XE-001
Output	Low-pressure gas discharge lines of xenon; diffuse aperture
Bulb life	~5000 hrs
Aperture	3.2 mm
Connector	905 SMA
Amplitude Stabilization	3 seconds
Nominal Frequency	28 kHz
Power Consumption	3 Watts
Power Requirements	5 VDC, 600 mA typical, 1500 mA max
Unit Dimensions	160 mm x 103 mm x 53 mm (6.30 in. x 4.06 in. x 2.09 in.)

Part Number: XE-001

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.