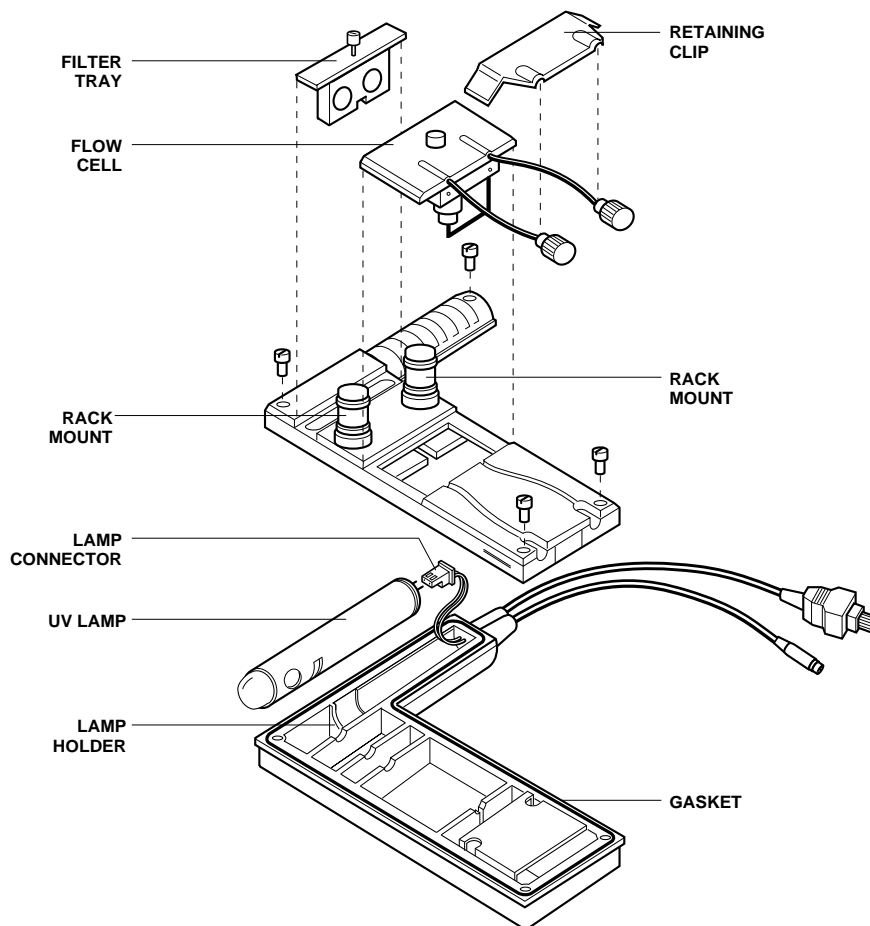


BioLogic LP Optics Module

Catalog Number 731-8324

When connected to the BioLogic LP system, the BioLogic LP Optics module detects biomolecules in a fluid stream. The portability of the optics module enables it to be positioned close to the outlet of a chromatography column, thus minimizing system dead volume and remixing of peaks.



UV Optics Module

Optics Module Installation

1. Plug the smaller round connector into the receptacle marked "UV Optics" on the rear panel of the BioLogic LP case.
2. Plug the larger rectangular connector into the receptacle marked "UV Lamp" on the rear panel of the BioLogic LP case.
3. Mount the module onto the rack using either one of the rack mounts located on the underside of the module.
4. Using male luer fittings, connect the inlet port of the optics module to the column outlet using the shortest amount of tubing possible. The direction of flow through the flow cell is indicated by raised arrows on top of the optics module.
5. Connect the exit port of the optics module to a fraction collector, diverter valve, or waste receptacle. If fractions are to be collected, it is best to minimize the diameter and length of the tubing used for connections.

Wavelength Selection

The BioLogic LP Optics Module has a filter tray containing both 254nm and 280nm filters (note markings on the filter unit). To change from one wavelength to another, loosen the thumbscrew and remove the filter tray. Rotate the filter tray 180° and reinsert, being careful not to over-tighten the thumbscrew. The selected wavelength is indicated by the raised arrow on the bottom of the case.

Flow Cell Replacement

WARNING: To avoid exposure to UV radiation, make sure that the UV lamp is off when changing filters and flow cells.

Flow cell specifications:

- 2mm path length
- Internal volume of 80 mL
- Illuminated volume of 3 µL

1. Remove the retaining clip from the bottom of the optics module and pull the luer connectors away from the case.
2. Unscrew the thumbscrew and remove the old flow cell.
3. Secure the o-ring on the replacement flow cell.
4. Reinsert the new flow cell following the above steps in the reverse order, being careful not to over-tighten the thumbscrew.

UV Lamp Replacement

1. Unplug the optics module from the rear panel of the BioLogic LP Controller.

WARNING: Failure to unplug the optics module from the control unit may expose the user to high voltage, UV radiation, and hot surfaces.

2. If the UV lamp has been on, allow the optics module to cool for ten minutes before continuing.
3. Remove the flow cell as described above. Remove the filter tray and the flow cell by unscrewing the thumbscrews then pulling the filter tray from the optics module.
4. Remove the four screws holding the optics module together and remove the bottom half of the case.
5. Pull the old UV lamp out of its holder and unplug it from its connector.

6. Using a protective cloth, plug the new UV lamp into its connector and insert it into the lamp holder.

NEVER handle the quartz surface of the UV lamp as grease and fingerprints will damage the lamp.

7. Reassemble the optics module, being careful that the gasket seal between the case halves is properly seated in its groove.

NOTE: An improperly seated gasket will result in light leaks, which will interfere with the proper operation of the instrument. One check for light leaks is to squeeze the two halves of the module and look for UV absorbance changes.

8. Replace the four screws holding the optics module together, being careful not to overtighten the screws.
9. Replace and secure the flow cell, filter tray, and retaining clip.

Ordering Information

731-8223	BioLogic LP Optics Module, replacement
731-8165	Flow Cell
731-8166	Lamp, Replacement
731-8167	Filter assembly, 280 & 254 nm
731-8155	Conductivity Flow Cell

BIO-RAD

Bio-Rad
Laboratories

Life Science
Group

Bio-Rad Laboratories Main Office, 2000 Alfred Nobel Drive, Hercules, California 94547, Ph. (510) 741-1000, Fx. (510) 741-5800
Also in: Reagents Park, Australia, Ph. 02-9914-2800, Fx. 02-9914-2888 Wien, Austria, Ph. (1) 877 89 01, Fx. (1) 876 56 29 Nazareth, Belgium, Ph. 09-385 55 11, Fx. 09-385 65 54
Mississauga, Canada, Ph. (905) 712-2771, Fx. (905) 712-2990 Beijing, China, Ph. (01) 2046622, Fx. (01) 2051876 Copenhagen, Denmark, Ph. 39 17 9947, Fx. 39 27 1698
Espoo, Finland, Ph. 90 804 2200, Fx. 90 804 1100 Ivry sur Seine Cedex, France, Ph. (1) 49 60 68 34, Fx. (1) 46 71 24 67 München, Germany, Ph. 089 31884-0, Fx. 089 31884-100
New Delhi, India, Ph. 91-11-461-0103, Fx. 91-11-461-0765 Milano, Italy, Ph. 02-21609.1, Fx. 02-21609.399 Tokyo, Japan, Ph. 03-5811-6270 Fx. 03-5811-6272 Veendam,
The Netherlands, Ph. 0318-540666, Fx. 0318-542216 Auckland, New Zealand, Ph. 09-443 3099, Fx. 09-443 3097 Kowloon, Hong Kong, Ph. 7893300, Fx. 7891257 Singapore,
Ph. (65) 272-9877, Fx. (65) 273-4835 Solna, Sweden, Ph. 46 (0) 8 735 83 00, Fx. 46 (0) 735 54 60 Madrid, Spain, Ph. (91) 661 70 85, Fx. (91) 661 96 98 Glatbrugg, Switzerland,
Ph. 01/809 55 55, Fx. 01/809 55 00 Hemel Hempstead, United Kingdom, Ph. 0800 181134, Fx. 01442 259118