

Bio-Rad Laboratories  
Improving lives with every test.

# Legionella

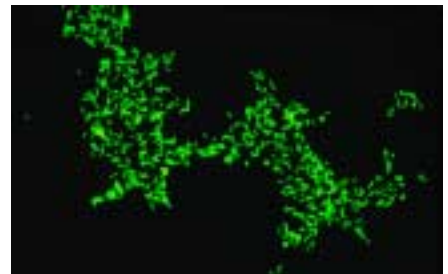
## MonoFluo<sup>®</sup> Legionella

**Several *Legionella* species are associated with human disease. More than 80% of *Legionella* organisms isolated from patients have been identified as *L. pneumophila*.**

*Legionella* are Gram-negative bacteria that are ubiquitous in the environment and may be present in fresh water lakes, tap water, air conditioning systems and recreational whirlpools.

Infections with *L. pneumophila* occur as sporadic, isolated events, endemic clusters of cases and epidemics. Transmission of the disease appears to be related to exposure to aerosols, such as those created by air conditioners, jet nebulizers and room humidifiers. It has been estimated that *L. pneumophila* accounts for approximately 10% of atypical pneumonias. Mortality rates for patients with Legionnaires' disease range from 15% of otherwise healthy individuals to 80% in untreated immunosuppressed individuals.

Clinical lab diagnosis to determine appropriate treatment is best done by direct examination of clinical specimens and by culture and isolation of *Legionella* from clinical specimens.



### Convenience

- Ready-to-use reagents and slides
- Monoclonal antibodies react to all known serogroups of *L. pneumophila*
- Rapid turnaround time

### Reliable

- 100% sensitivity
- 100% specificity

*MonoFluo Legionella pneumophila IFA Test Kit*

**BIO-RAD**



# Legionella Testing

## Immunofluorescence

The MonoFluo® *Legionella pneumophila* IFA Test kit is a user-friendly indirect immunofluorescence assay with ready-to-use reagents in dropper bottles. A single monoclonal antibody is used to detect and identify all serogroups of *L. pneumophila* directly in patient specimens or in culture isolates, confirming a diagnosis of Legionnaires' disease.

## MonoFluo® *Legionella pneumophila* IFA Test Kit

Component	Contents	Preparation
R1 • Anti- <i>Legionella pneumophila</i> Staining Reagent 1 dropper bottle (1.3 mL)	<ul style="list-style-type: none"> <li>FITC-Labeled Monoclonal Antibody</li> <li>Counterstain</li> <li>Sodium Azide</li> <li>Protein-stabilized buffer</li> </ul>	Use as supplied
C1 • Positive Control Antigen Suspension 1 dropper bottle (2.0 mL)	<ul style="list-style-type: none"> <li>Killed <i>L. pneumophila</i></li> <li>Buffered Saline</li> <li>Sodium Azide</li> </ul>	Shake before use
R2 • Mounting Medium 1 dropper bottle (4.0 mL)	<ul style="list-style-type: none"> <li>Buffered glycerol</li> <li>Anti-Quencher</li> </ul>	Use as supplied
R3 • Fluorescence Microscopy Slides 24 (2 wells each)	<ul style="list-style-type: none"> <li>Fluorescence microscopy slides</li> </ul>	Wipe with lint-free tissue before use
Store reagents at 2-8°C. Slides should be stored at 2-28°C.		

## Ordering Information

Catalog No.	Description
32514	MonoFluo® <i>Legionella pneumophila</i> IFA Test Kit.....24 tests



Bio-Rad Laboratories

For further information, please contact the Bio-Rad office nearest you or visit our website at [www.bio-rad.com](http://www.bio-rad.com).

Clinical  
Diagnostics Group

Website [www.bio-rad.com](http://www.bio-rad.com) U.S. 1-800-2-BIO-RAD Australia 61-2-9914-2800 Austria 43-1-877-8901 Belgium 32-9-385-5511 Brazil 5521-2507-6191 Canada 1-514-334-4372 Czech Republic 420-2-41430532 China 86-21-63052255 Denmark 45-4452-1000 Finland 358-9-804-22-00 France 33-1-4795-6000 Germany 49-89-31884-0 Hong Kong 852-2789-3300 India 91-124-6398112 Israel 972-3-9514127 Italy 39-02-216091 Japan 81-3-5811-6290 Korea 82-2-3473-4460 Latin America 305-894-5950 Mexico 5255-5534-2552 The Netherlands 31-318-540666 New Zealand 64-9-415-2280 Norway 47-23-38-41-30 Poland 48-22-331-99-99 Portugal 351-21-472-7700 Russia 7-095-721-14-00 Singapore 65-6415-3188 South Africa 27-11-442-85-08 Spain 34-91-590-5200 Sweden 46-8-555-127-00 Switzerland 41-61-717-95-55 Thailand 662-651-8311 United Kingdom 44-208-328-2000