



# IT LEISH Individual Rapid Test

Leishmaniasis Diagnosis, Anywhere



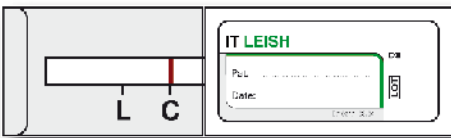
# IT LEISH Individual Rapid Test a Reference Tool for Visceral Leishmaniasis Diagnosis

Leishmaniasis is a vector-borne disease transmitted by phlebotomine sandflies in many tropical and sub-tropical countries. Visceral leishmaniasis (VL) is the most serious form of the disease, and has an estimated incidence of 500.000 new cases and 60.000 deaths per year. Most cases of VL are found in north-eastern India area, eastern Africa and South-America. VL is fatal if left untreated, with a fatality rate in developing countries that can reach 100% within 2 years. Early diagnosis is considered an essential component of VL control. VL diagnosis tests used in the field are based on anti-*Leishmania* antibodies detection. Among them, immunochromatographic rapid tests based on rk39 protein are mostly used.

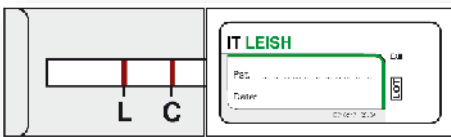
- **IT LEISH** is a reference immuno-chromatographic RDT using **rk39 antigen** to confer high specificity and high sensitivity of detection.
- **IT LEISH** is an accurate assay for **confirmation of clinically suspected cases** and early **initiation of adequate treatment**.
- **IT LEISH** is individually packed and includes all components required for whole blood sampling and test processing. The test can be performed in **any field situation** and after a **short educational training period**.
- **IT LEISH** results are obtained in **20 minutes** only.

## Result Interpretation (see Product Insert)

### NEGATIVE



### POSITIVE



## Bibliography

- Evaluation of a new recombinant k39 rapid diagnostic test for Sudanese visceral leishmaniasis. Ritmeijer K. and *al*, Am. J. Trop. Med. Hyg., 2006.
- Rapid, noninvasive diagnosis of visceral leishmaniasis in India: comparison of two immunochromatographic strip tests for detection of anti-K39 antibody. Sundar S. and *al*, J. Clin. Microbio., 2006.
- Field evaluation of rk39 test and direct agglutination test for diagnosis of visceral leishmaniasis in a population with high prevalence of human immunodeficiency virus in Ethiopia. Horst R. and *al*, Am. J. Trop. Med. Hyg., 2009.

## Ordering Information

Catalog No.	Description
710124	IT LEISH Individual Rapid Test ..... 24 tests/kit



Bio-Rad  
Laboratories

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

Clinical  
Diagnostics Group

Website [www.bio-rad.com/diagnostics](http://www.bio-rad.com/diagnostics) U.S. 1-800-2BI/O-RAD Australia 61-2-9914-2800 Austria 43-1-877-8901 Belgium 32-9-385-5511 Brazil 5521-3237-9400 Canada 1-514-334-4372 China 86-21-64260808 Czech Republic 420-241-430-532 Denmark +45-4452-1000 Finland 358-9-804-22-00 France 33-1-47-95-60-00 Germany +49-(0)89-318-840 Greece 30-210-7774396 Hong Kong 852-2789-3300 Hungary +36-1-459-6100 India 1-800-180-1224 Israel 972-3-9636050 Italy +39-02-216091 Japan 81-3-6361-7070 Korea 82-2-3473-4460 Mexico +52 (55) 5488-7670 The Netherlands +31-318-540666 New Zealand 64-9-415-2280 Norway 47-23-38-41-30 Poland 48-22-3319999 Portugal 351-21-472-7700 Russia 7-495-721-14-04 Singapore 65-6415-3170 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-(0)20-8328-2000