



QIAstat-Dx Meningitis/Encephalitis Panel

For quick and accurate detection of central nervous system infections



Meningitis and encephalitis are medical emergencies. Rapid diagnosis and treatment initiation are crucial for successful patient outcomes. Traditional diagnostic methods alone involve multiple time-consuming tests and can take 2 days or more. Imagine getting 15 results in a little over an hour. Powered by our easy workflow and reliable multiplex PCR technology, the QIAstat-Dx Meningitis/Encephalitis Panel makes this possible.

- **Get comprehensive results fast**

- Detect and differentiate 15 pathogen targets in about an hour
- Support timely clinical decision-making and rapid therapeutic intervention

- **Set up your tests quickly and easily**

- All-in-one assay cartridges require less than one minute hands-on time
- Just 200 μ L cerebrospinal fluid with no precision pipetting required

- **Receive supportive information**

- Easily access Ct values and amplification curves for all detected pathogens



Comprehensive pathogen coverage with QIAstat-Dx

Bacterial

- *Escherichia coli* K1
- *Haemophilus influenzae*
- *Listeria monocytogenes*
- *Neisseria meningitidis* (encapsulated)

- *Streptococcus agalactiae*
- *Streptococcus pneumoniae*
- *Mycoplasma pneumoniae*
- *Streptococcus pyogenes*

Viral

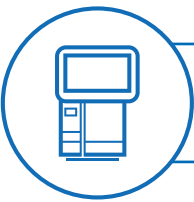
- Enterovirus
- Herpes simplex virus 1
- Herpes simplex virus 2
- Human herpes virus 6
- Human parechovirus
- Varicella zoster virus

Fungal

- *Cryptococcus neoformans/gattii**

**Cryptococcus neoformans*/*Cryptococcus gattii* are both detected, but not differentiated.

High overall clinical sensitivity (95.24% (CI, 90.50–97.67%)) and specificity (99.93% (CI 99.84–99.97%))†



QIAstat-Dx Meningitis/Encephalitis Panel runs on our modular QIAstat-Dx Analyzer.

Ordering Information - Assay

Product	Contents	Cat. no.
QIAstat-Dx Meningitis/Encephalitis Panel	Six individually packaged cartridges containing all reagents needed for sample preparation and multiplex RT-real time PCR plus internal control, including six transfer pipettes	691611



Confidently detect central nervous system infections with QIAstat-Dx – learn more at [QIAGEN.com/QIAstat-Dx](https://www.qiagen.com/QIAstat-Dx)

† Data presented is for the QIAstat-Dx Meningitis/Encephalitis Panel (Cat. No. 691611); QIAstat-Dx Meningitis/Encephalitis Panel Instructions for Use. QIAGEN, April 2022

The QIAstat-Dx Analyzer is intended for in vitro diagnostic use.

The QIAstat-Dx ME Panel is indicated as an aid in the diagnosis of specific agents of meningitis and/or encephalitis and results must be used in conjunction with other clinical, epidemiological, and laboratory data. Results from the QIAstat-Dx ME Panel are not intended to be used as the sole basis for diagnosis, treatment, or other patient management decisions. Positive results do not rule out co-infection with organisms not included in the QIAstat-Dx ME Panel. The agent or agents detected may not be the definite cause of the disease. Negative results do not preclude central nervous system (CNS) infection. The QIAstat-Dx ME Panel is intended for in vitro diagnostic use by laboratory professionals only.

Trademarks: QIAGEN®, Sample to Insight®, QIAstat-Dx® (QIAGEN Group). PROM-15028-002 1128510 07/2022 © 2022 QIAGEN, all rights reserved.

Ordering
Technical Support
Website

www.qiagen.com/shop
www.support.qiagen.com
www.qiagen.com