



## QIAstat-Dx Gastrointestinal Panel 2

# For quick and accurate detection of gastrointestinal infections



Traditional diagnostic methods for gastrointestinal infections, including bacterial culture, can be slow and cumbersome. They often aren't very sensitive and provide low diagnostic yield. Imagine getting 22 results in about an hour. Powered by our easy workflow and reliable multiplex PCR technology, the QIAstat-Dx Gastrointestinal Panel 2 makes this possible.

- **Get comprehensive results fast**

- Detect and differentiate 22 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  $\mu$ L liquid transport medium 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

- *Campylobacter* (*C. jejuni*, *C. upsaliensis*, *C. coli*)
- *Clostridium difficile* (toxin A/B)
- Enteroaggregative *E. coli* (EAEC)
- Enteroinvasive *E. coli* (EIEC)/*Shigella*
- Enteropathogenic *E. coli* (EPEC)
- Enterotoxigenic *E. coli* (ETEC) *lt/st*
- *Plesiomonas shigelloides*
- *Salmonella* spp.
- Shiga-like toxin-producing *E. coli* (STEC) *stx1/stx2*\* (including specific identification of *E. coli* O157 sero-group within STEC)
- *Vibrio vulnificus*
- *Vibrio parahaemolyticus*
- *Vibrio cholerae*
- *Yersinia enterocolitica*

## Viral

- Adenovirus F40/41
- Astrovirus
- Norovirus (GI, GII)
- Rotavirus A
- Sapovirus (GI, GII, GIV, GV)

## Parasitic

- *Cryptosporidium*
- *Cyclospora cayetanensis*
- *Entamoeba histolytica*
- *Giardia lamblia*

\*Toxin content of STEC (*stx1* and/or *stx2*) is reported by the QIAstat-Dx Gastrointestinal Panel 2



QIAstat-Dx Gastrointestinal Panel 2 runs on both our modular **QIAstat-Dx Analyzer** and higher-throughput **QIAstat-Dx Rise**.

## Ordering Information - Assay

Product	Contents	Cat. no.
QIAstat-Dx Gastrointestinal Panel 2	Six individually packaged cartridges containing all reagents needed for sample preparation and multiplex RT-real time PCR plus internal control, including six transfer pipettes	691412



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

The QIAstat-Dx Analyzer, QIAstat-Dx Rise and QIAstat-Dx Gastrointestinal Panel 2 are intended for in vitro diagnostic use.

Trademarks: QIAGEN®, Sample to Insight®, QIAstat-Dx® (QIAGEN Group). PROM-20910-001 1127824 06/2022 © 2022 QIAGEN, all rights reserved.

Ordering  
Technical Support  
Website

[www.qiagen.com/shop](https://www.qiagen.com/shop)  
[www.support.qiagen.com](https://www.support.qiagen.com)  
[www.qiagen.com](https://www.qiagen.com)