



# Molecular Controls for STI Testing

**Reduce risk by increasing compliance.**

AmpliTrol™, AmpliPROBE™, and AmpliClear™ quality control products are designed to monitor the performance of nucleic acid testing of Sexually Transmitted Infections (STI). This family of products includes quality controls for *Chlamydia trachomatis* (CT), *Neisseria gonorrhoeas* (NG or GC), and Human Papillomavirus (HPV) assays.

**Guidelines recommend independent quality controls to minimize medically important analytical errors.**

Guidelines call for independent controls because instrument manufacturer controls are less sensitive to slight changes in device performance and could fail to detect errors that can potentially be harmful to patients.

*“Use of independent third party control materials should be considered, either instead of, or in addition to, any control materials supplied by the reagent or instrument manufacturer.”*

**ISO 15189, Medical Laboratories – Requirements for Quality and Competence**

## AmpliTrol™ CT/GC

For use as an unassayed quality control in DNA amplification assays detecting *Chlamydia trachomatis* and *Neisseria gonorrhoeae*.

- 2 year shelf life at 2–8°C
- 60 day open-vial stability at 2–8°C

### Analytes

*Chlamydia trachomatis*  
*Neisseria gonorrhoeae*

## AmpliPROBE™ CT/GC

For use as an unassayed quality control in the Aptima Combo 2® test system for detecting *Chlamydia trachomatis* and *Neisseria gonorrhoeae*.

- 2 year shelf life at 2–8°C
- 60 day open-vial stability at 2–8°C

### Analytes

*Chlamydia trachomatis*  
*Neisseria gonorrhoeae*

Refer to [myinserts.com](http://myinserts.com) or the package insert of currently available lots for specific analyte and stability claims. Please note that not all product may be available in all countries.



## AmpliClear™

For use as an unassayed, non-reactive quality control in nucleic acid amplification assays detecting *Chlamydia trachomatis* and *Neisseria gonorrhoeae*.

- 2 year shelf life at 2–8°C
- 60 day open-vial stability at 2–8°C

### Analytes

#### Non-reactive for:

*Chlamydia trachomatis*  
*Neisseria gonorrhoeae*

Refer to [myinserts.com](http://myinserts.com) or the package insert of currently available lots for specific analyte and stability claims.

## AmpliTrol™ HPV

For use as an unassayed quality control for detecting Human Papillomavirus.

- 2 year shelf life at 2–8°C
- 60 day open-vial stability at 2–8°C

### Analyte

Human Papillomavirus (HPV)

## Use Bio-Rad Molecular Quality Controls on these popular assays

- Available for Assay
  - Positive on type 16 HPV only
- Manufacturer/Assay(s)

	AmpliClear	AmpliPROBE CT/GC	AmpliTrol CT/GC	AmpliTrol HPV
<b>Abbott Diagnostics</b>				
m2000 RealTime CT/NG	●		●	
m2000 RealTime HPV				○
<b>Becton Dickinson</b>				
BD ProbeTec™ ET CT/NG	●		●	
<b>Cepheid®</b>				
Xpert® CT/NG			●	
<b>Hologic®</b>				
APTIMA® Combo 2®	●	●		
APTIMA® HPV				○
<b>QIAGEN</b>				
digene HC2 HPV DNA Test				●
<b>Roche Diagnostics</b>				
cobas® 4800 CT/NG	●		●	
cobas® 4800 HPV				○

If your assay is not listed, please contact your nearest Bio-Rad office for current information.

## Ordering Information

<b>Cat #</b>	<b>Configuration</b>	<b>Cat #</b>	<b>Configuration</b>
<b>AmpliTrol™ CT/GC</b>		<b>AmpliClear™</b>	
00126	Single Level/Dropper Bottle . . . . . 1 x 5 mL	00127	Single Level/Dropper Bottle . . . . . 1 x 5 mL
<b>AmpliPROBE™ CT/GC</b>		<b>AmpliTrol™ HPV</b>	
00138	Single Level/Dropper Bottle . . . . . 1 x 5 mL	00139	Single Level/Dropper Bottle . . . . . 1 x 5 mL



**Bio-Rad Laboratories**

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

### Clinical Diagnostics Group

Website [www.bio-rad.com/qualitycontrol](http://www.bio-rad.com/qualitycontrol) Australia 61-2-9914-2800 Austria 43-1-877-8901 Belgium +32 (3)710-53-00 Brazil +55 (31)3689-6600 Canada 1-514-334-4372 China 86-21-61698500 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 1800-180-1224 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 +7-495-721-1404 Singapore 65-6415-3170 South Africa 27-11-442-85-08 Spain 34-91-590-5200 Sweden +46-8-555-127-00 Switzerland +41 (0)26-674-55-05/06 Taiwan 886-2-2578-7189 Thailand 662-651-8311 United Kingdom +44 (0)20-8328-2000

ProbeTec is a trademark of Beckton, Dickinson and Company. Cepheid and Xpert are registered trademarks of Cepheid. APTIMA and Combo 2 and Hologic are trademarks of Hologic. COBAS is a trademark of Roche Molecular Systems Inc.