



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 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 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
Abbott Diagnostics				
m2000 RealTime CT/NG	●		●	
m2000 RealTime HPV				○
Becton Dickinson				
BD ProbeTec™ ET CT/NG	●		●	
Cepheid®				
Xpert® CT/NG			●	
Hologic®				
APTIMA® Combo 2®	●	●		
APTIMA® HPV				○
QIAGEN				
digene HC2 HPV DNA Test				●
Roche Diagnostics				
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

Cat #	Configuration	Cat #	Configuration
AmpliTrol™ CT/GC		AmpliClear™	
00126	Single Level/Dropper Bottle 1 x 5 mL	00127	Single Level/Dropper Bottle 1 x 5 mL
AmpliPROBE™ CT/GC		AmpliTrol™ HPV	
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**

**Clinical
Diagnostics Group**

Website 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.