

Immunoassay Controls

Discover a full range of quality control products for your reproductive and prenatal testing needs





Bio-Rad offers a comprehensive quality control solution for today's reproductive health and prenatal testing

An estimated 48.5 million couples worldwide are unable to conceive a child after five years of trying to become pregnant.¹ In the United States, the quality of reproductive health testing has never been more important. The Center for Disease Control reports that an estimated 7.3 million women use fertility services, indicating that infertility is a developing concern.² Bio-Rad can help alleviate that concern by providing reliable monitoring of fertility testing.

Bio-Rad quality controls cover a comprehensive menu of analytes designed to fit your reproductive health and fertility testing needs. For greater confidence in reliability of patient test results, manage your QC data and access peer group comparisons with our Unity Interlaboratory Program.

¹ Mascarenhas, M.N., Flaxman, S.R., Boerma, T., Vanderpoel, S., Stevens, G.A., "National, Regional, and Global Trends in Infertility Prevalence Since 1990: A Systematic Analysis of 277 Health Surveys." *PLoS Med* 9(12): e1001356 (2012).

² Centers for Disease Control and Prevention (CDC), National Center for Health Statistics, Key Statistics from the National Survey of Family Growth, (data 2011-2015)



Reproductive Health Quality Controls

Liquichek AMH Control



A liquid, assayed quality control serum with clinically significant levels for monitoring ovarian reserve test procedures. Choose an ideal quality control solution for your clinic or laboratory.

- Liquid convenience to reduce handling errors
- Three levels targeted at low, medium and high clinically-relevant ranges
- Access to the Unity Interlaboratory Program, powerful and reliable quality control data management
- Human base to more closely resemble a patient sample
- 26 month shelf life at -20°C to -70°C
- 14 day open-vial stability at 2-8°C

Analyte

Anti-Mullerian Hormone

Lyphochek Fertility Control



- Designed to monitor the precision of tests performed for fertility or pregnancy
- Elevated hCG and hCG-β Subunit levels allow laboratories to monitor their dilution procedure
- Human serum based
- 3 year shelf life at 2-8°C
- 7 day reconstituted stability at 2-8°C

POCT Methods Listed

Analytes

Alphafetoprotein (AFP)
Digitoxin
Estradiol
Estril (Free)
Estril (Total)*
Ferritin
FSH

hCG
hCG-β Subunit
LH
Progesterone
Testosterone

*No claims are made regarding performance or stability.

Liquichek Immunoassay Plus Control



- Offers a vast array of popular routine immunoassay analytes
- Assigned values are available for all major automated analyzers
- Provides a highly efficient solution for laboratories that focus on routine tests
- Liquid, human serum based
- 2 year shelf life at -20°C to -70°C
- 14 day open-vial stability at 2-8°C for most analytes

POCT Methods Listed

Analytes

For Reproductive Health Testing

Androstenedione
DHEA Sulfate
Estradiol
FSH
LH
Progesterone
Prolactin
SHBG (Sex Hormone Binding Globulin)
Testosterone

Prenatal Testing Quality Controls

Liquichek Maternal Serum II Control



- Liquid, assayed quality control serum
- Analytes at clinically significant levels to monitor the precision of maternal serum analysis
- Human serum
- Includes the hormone Inhibin A
- Very elevated hCG
- Very low levels of Free Estriol and AFP

Analytes

AFP
hCG
Free Estriol
Inhibin A

Liquichek Maternal Serum 1st Trimester Control



- Liquid, assayed quality control
- Designed to monitor the precision of immunoassay test methods used during maternal serum first trimester screening
- Includes clinically relevant levels of Free β -hCG and PAPP-A
- Human serum based
- 2 year shelf life at -20°C to -70°C
- 15 day open-vial stability at $2-8^{\circ}\text{C}$

Analytes

hCG
hCG- β Subunit (Free)
PAPP-A

Available outside the United States only.

Essential tests for reproductive health available in Bio-Rad quality controls

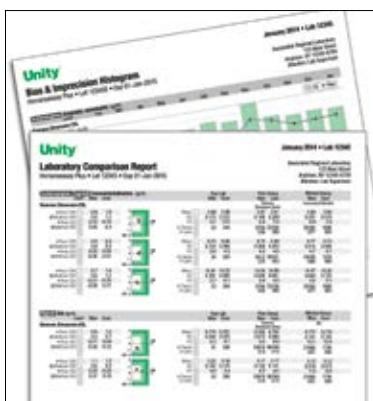
Test	Evaluation of Infertility		Liquichek Immunoassay Plus Control	Liquichek AMH Control	Lyphocheck Fertility Control
	Women	Men			
AMH	●			●	
Androstenedione	●	●	●		
DHEAS	●	●	●		
Estradiol	●	●	●		●
FSH	●	●	●		●
LH	●		●		●
Progesterone	●		●		●
Prolactin	●	●	●		
SHBG	●	●	●		
Testosterone	●	●	●		●
TSH	●		●		



Unity QC Data Management Solutions

Bio-Rad Quality Controls are part of the Unity Program with its industry-leading features for benchmarking laboratory performance.

- Benefit from peer data generated from more than 50,000 instruments worldwide
- Utilize real-time access to peer group comparisons to help troubleshooting
- Consolidate your entire laboratory into one system
- Facilitate run-time decisions with optimized Westgard rules
- Ease compliance with ISO 15189 using simple tools to review rule violations and document corrective actions
- Define your own quality requirements and track key quality performance indicators



You can consolidate your entire laboratory into one QC system through our Unity Program. This sophisticated software includes best-in-class tools to support releasing results, benchmarking laboratory performance, and designing your QC system.

Unity Software Solutions

Contact your local sales representative for more information about our range of QC Data Management Solutions



Unity Real Time

Expert QC Data Management Solution for Desktop or Online Users

- Facilitate CLIA and ISO 15189 compliance
- Improve real-time bench and supervisor QC data review
- Implement best QC rules when used with Westgard Advisor
- Run validation with comprehensive audit trails
- Advanced charts and reports for data analysis
- Reduce non-essential retests with Analytical Goal options
- Upload QC data points from an LIS, middleware or instrument (optional)
- RiLiBÄK Advisor module available to comply with German regulations

Ordering Information

Cat # Description

Liquichek AMH Control

12006327 Trilevel (2 of each level) 6 x 2 mL

Lyphochek Fertility Control

480 Trilevel (4 of each level) 12 x 5 mL
 481 Level 1 12 x 5 mL
 482 Level 2 12 x 5 mL
 483 Level 3 12 x 5 mL
 480X Trilevel MiniPak (1 of each level)3 x 5 mL

Liquichek Immunoassay Plus Control

360 Trilevel (4 of each level) 12 x 5 mL
 361 Level 1 12 x 5 mL
 362 Level 2 12 x 5 mL
 363 Level 3 12 x 5 mL
 360X Trilevel MiniPak (1 of each level)3 x 5 mL

Liquichek Maternal Serum II Control

402 Level 1 6 x 2.5 mL
 403 Level 2 6 x 2.5 mL
 404 Level 3 6 x 2.5 mL
 403X Trilevel MiniPak (1 of each level) 3 x 2.5 mL

Liquichek Maternal Serum 1st Trimester Control*

636 Trilevel (2 of each level)6 x 1 mL
 636X Trilevel MiniPak (1 of each level)3 x 1 mL

*Available outside the US only

Unity Real Time

804-1 Unity Real Time. Annual subscription

Bio-Rad Mission: Control

1100002 Bio-Rad Mission: Control Annual Subscription
 1100005 Bio-Rad Mission: Control Subscription & Service Bundle



EQAS

An independent, external assessment of performance in comparison to your peers.



Independent QC

Ongoing, proactive, unbiased daily QC that helps identify errors as they occur or begin to trend.



Unity

QC Data Management tools that help you create a strategy to reduce risk and streamline QC workflow.



**Bio-Rad
Laboratories, Inc.**

For further information, please contact the Bio-Rad office nearest you or visit our website at www.bio-rad.com/iacontrols

Clinical
Diagnostics Group

Website www.bio-rad.com/qualitycontrol USA 1 800 224 6723 Australia 61 2 9914 2800 Austria 43 01 877 89019 Belgium 32 03 710 53 00 Brazil 55 11 3065 7550
 Canada 1 905 364 3435 China 86 21 6169 8500 Czech Republic 36 01 459 6192 Denmark 45 04 452 10 00 Finland 35 08 980 422 00 France 33 01 479 593 00
 Germany 49 089 3188 4393 Hong Kong 852 2789 3300 Hungary 36 01 459 6190 India 91 124 4029300 Israel 972 03 963 6050 Italy 39 02 49486600 Japan 81 3 6361 7000
 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 310 318 540 666 New Zealand 64 9 415 2280 Norway 47 0 233 841 30 Poland 36 01 459 6191
 Portugal 351 21 4727717 Russia 7 495 721 14 04 Singapore 65 6415 3188 South Africa 36 01 459 6193 Spain 34 091 49 06 580 Sweden 46 08 555 127 00
 Switzerland 41 0617 17 9555 Taiwan 886 2 2578 7189 Thailand 66 2 651 8311 United Arab Emirates 971 4 8187300 United Kingdom 44 01923 47 1301

Bio-Rad, Liquichek, Lyphochek and Unity Real Time are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions.

© 2018 Bio-Rad Laboratories, Inc. Printed in USA QSD17-098 Q-1659 Rev. 04/2018

Printed on recycled paper with soy-based inks