

Liquichek™ Maternal Serum 1st Trimester Control

A liquid, trilevel control designed to monitor the precision of maternal serum analysis. Includes analytes commonly tested during first trimester screening.

- Human serum based
- 2 year shelf life at -20°C to -70°C
- 15 day open-vial stability at 2–8°C



Analytes

hCG
hCG-β Subunit (Free)
PAPP-A

Refer to the package insert of currently available lots for specific analyte and stability claims

Ordering Information

Cat #	Liquichek™ Maternal Serum 1st Trimester Control	Quantity
636	Trilevel (2 of each level)	6 x 1 mL
636X	Trilevel MiniPak (1 of each level)	3 x 1 mL

Available outside the United States only.

Bio-Rad is Dedicated to Advancing QC Solutions Worldwide

Quality Controls

Broad portfolio of independent controls provides an unbiased assessment of methods and system performance.

QC Data Management Solutions

Manage and interpret your QC results using the industry-leading Unity™ Interlaboratory Program and benefit from data generated by users worldwide.

Customer Support

Receive the backing of full-service technical and customer support trained to handle your inquiries in a prompt, knowledgeable manner.

Education

We offer the educational tools and resources to help improve your ability to produce reliable test results.

Efficiently manage your QC product inserts with My eInserts™. Visit www.myeinserts.com for more information.



**Bio-Rad
Laboratories**

*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 **Australia** 61-2-9914-2800 **Austria** 43-1-877-8901 **Belgium** 32-9-385-5511 **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** 7-495-721-14-04 **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