



Liquichek™ Maternal Serum II Control



Liquichek™ Maternal Serum II Control

A liquid, assayed quality control serum with 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

Refer to myinserts.com or the package insert of currently available lots for specific analyte and stability claims.

Ordering Information

Cat #	Liquichek™ Maternal Serum II Control	Configuration
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



**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 (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