



# Liquichek™ Tumor Marker Control



## Liquichek™ Tumor Marker Control

Liquichek™ Tumor Marker Control is a liquid, human serum based, third party control for monitoring the precision of tumor marker testing procedures. This multi-analyte control has 3 levels and features 17 cancer antigens and popular tumor markers. The Unity™ Interlaboratory Program with peer group comparison is available for use with this control.



- Liquid convenience, human serum based, three desirable levels
- Third party control provides an unbiased assessment of the test systems
- Multi-analyte (17 tumor markers)
- Very high levels of cancer antigens (CA) and Ferritin
- Very low PSA level for highly sensitive PSA test methods
- Excellent 30 day open-vial stability at 2–8°C for most analytes
- 2 year shelf life when stored frozen (-20 to -70°C); minimizes frequency of crossover studies
- Assayed values available for most manual and automated test methods
- Peer group reports available through Unity™ Interlaboratory Program

### Analytes

AFP	CA 27.29	Cyfra 21-1*	PAP	PSA (Free)
2-Microglobulin	CA 72-4*	Ferritin	Prolactin	Thyroglobulin
CA 15-3	CA 125	hCG/ -hCG	PSA	
CA 19-9	CEA	IGF-1		

\*International Use Only. Not available for diagnostic use in the United States.

### Ordering Information

Cat #	Liquichek™ Tumor Marker Control	Quantity
547	Level 1 .....	6 x 2 mL
548	Level 2 .....	6 x 2 mL
549	Level 3 .....	6 x 2 mL
548X	Trilevel MiniPak (1 of each level).....	3 x 2 mL



This control is featured in the Unity™ Interlaboratory Program. Participation in the Unity™ Interlaboratory Program can help improve your laboratory quality and help you meet regulatory and accreditation requirements worldwide. You may also subscribe to one of our Unity™ QC Data Management Solutions designed to improve the effectiveness of your statistical process control. Ask your Bio-Rad Account Representative for more information.

**Discover the power of Unity™ Solutions at [www.QCNet.com](http://www.QCNet.com).**



**Bio-Rad  
Laboratories**

*For further information, please contact your local Bio-Rad office  
or learn more online at [www.bio-rad.com/qualitycontrol](http://www.bio-rad.com/qualitycontrol)*

*Clinical  
Diagnostics Group*

**Website** [www.bio-rad.com/qualitycontrol](http://www.bio-rad.com/qualitycontrol) **U.S.** 1-800-2BIO-RAD **Australia** 61-2-9914-2800 **Austria** 43-1-877-8901 **Belgium** 32-9-385-5511 **Brazil** 5521-3237-9400  
**Canada** 1-514-334-4372 **China** 86-21-64260808 **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** 36-1-455-8800 **India** 91-124-4029300 **Israel** 972-3-9636050  
**Italy** +39-02-216091 **Japan** 81-3-6361-6902 **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** 65-6415-3188 **South Africa** 27-11-442-85-08  
**Spain** 34-91-590-5200 **Sweden** 46-8-555-127-00 **Switzerland** 41-61-717-95-55 **Thailand** 662-651-8311 **United Kingdom** +44-(0)20-8328-2000