



Cardiac Assessment Controls with Values for Roche Instrument Systems

Introducing Liquichek™ Cardiac Troponins Control
with values for Troponin T hs (outside the US only)





Quality Control for Troponin I and Troponin T

Values for Troponin T hs (high sensitive)*

Liquichek™ Cardiac Troponins Control

Quality Control of Troponin assays at low levels is critical because results above the 99th percentile upper reference limit (URL) designate myocardial infarction.** Liquichek™ Cardiac Troponins Control has been developed to meet the urgent demand from Roche customers for control of Troponin T hs at this key medical decision point.

Roche cobas® Assay	Medical Decision Point	Liquichek™ Cardiac Troponins Control	Roche PreciControl Troponin
Troponin I	0.3 ng/mL	0.5	0.75
Troponin T hs	14 pg/mL	14	30

International Organization for Standardization (ISO) 15189 recommends the use of independent control materials either instead of, or in addition to, any control materials supplied by the reagent or instrument manufacturers.

Liquichek™ Cardiac Troponins Control compared to Roche PreciControl Troponin:

- Values closer to the Medical Decision Point
- Convenience of a liquid control
- Significantly longer open-vial stability (longer shelf life coming soon)

Analytes

Troponin I
Troponin T

Control Comparison	Liquichek™ Cardiac Troponins Control	Roche PreciControl Troponin
Form	Liquid	Lyophilized
Open-vial stability for Troponin T	25 days at 2-8°C	4 days at 2-8°C
Shelf life	18 months at -20 to -70°C	~18 months at 2-8°C
Packaging	6 x 3 mL (MiniPak also available)	4 x 2 mL (2 of each level)

For higher values of Troponin I and Troponin T, use Levels 2 and 3 of Liquichek™ Cardiac Markers Plus Control LT.

*Values for Troponin T hs are provided for use outside U.S. only. U.S. customers are provided values for Troponin I and Troponin T (Gen 4).

**Joint ESC/AACF/AHA/WHF Task Force for the Universal Definition of Myocardial Infarction. Third universal definition of myocardial infarction. European Heart Journal 2012; 33: 2551-2567. Available online at: www.escardio.org/guidelines.

For specific analyte, open-vial stability, methods, mean values and ranges, refer to the package inserts of currently available lots at myeinserts.com



Quality Control for additional Cardiac Markers

Liquichek™ Cardiac Markers Plus Control LT

A four level cardiac assessment control designed to monitor assay precision across a wide range of values.

- Liquid, human serum based
- 3 year shelf life at -20°C to -70°C
- 20 day open-vial stability for most analytes
- Values provided for Troponin T (Gen 4)
- Level Low offers lower target values for all analytes

Analytes

BNP
 CK (Total)
 CK-MB Isoenzyme
 Digitoxin
 Homocysteine*
 hs-CRP
 Myoglobin
 NT-proBNP
 Troponin I
 Troponin T

*No claim is made for expected values or stability.

Liquichek™ D-dimer Control

A four level, single analyte control designed to monitor D-dimer assay precision across a wide assay range.

- Liquid, human plasma based
- 3 year shelf life at -20°C to -70°C
- 15 day open-vial stability at 2-8°C

Analyte

D-dimer

Liquichek™ Homocysteine Control

A trilevel, single analyte control designed to monitor Homocysteine precision across a wide assay range.

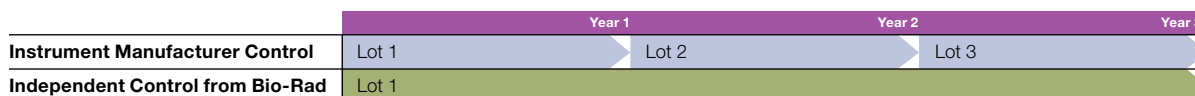
- Liquid, human serum based
- 3 year shelf life at -20°C to -70°C
- 30 day open-vial stability at 2-8°C

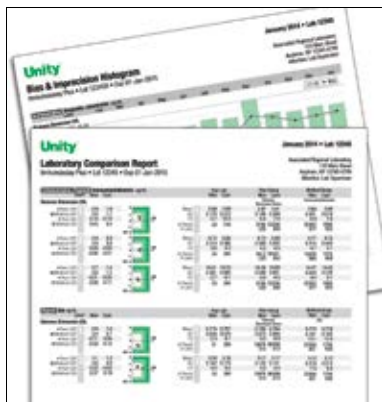
Analyte

Homocysteine

Save time and money through lot reservation

Perform fewer cross-over studies by reserving lots for up to three years.





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 40,000 instruments worldwide
- Utilize real-time access to peer group comparisons to help with troubleshooting
- Consolidate your entire laboratory into one QC system
- Facilitate run-time decisions with optimized Westgard rules
- Facilitate compliance with 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.

Ordering Information

Cat #	Product Name	Quantity	Cat #	Product Name	Quantity
Liquichek™ Cardiac Troponins Control			Liquichek™ D-dimer Control		
12001084	Single Level	6 x 3 mL	27100	Level Low	6 x 1 mL
12001085	MiniPak	1 x 3 mL	27101	Level 1	6 x 1 mL
Liquichek™ Cardiac Markers Plus Control LT			27102	Level 2	6 x 1 mL
146	Level 1	6 x 3 mL	27103	Level 3	6 x 1 mL
147	Level 2	6 x 3 mL	27102X	Four Level MiniPak (1 of each level)	4 x 1 mL
148	Level 3	6 x 3 mL	Liquichek™ Homocysteine Control		
149	Level Low	6 x 3 mL	287	Level 1	6 x 1 mL
146X	Level 1 MiniPak	1 x 3 mL	288	Level 2	6 x 1 mL
147X	Level 2 MiniPak	1 x 3 mL	289	Level 3	6 x 1 mL
148X	Level 3 MiniPak	1 x 3 mL	288X	Trilevel MiniPak (1 of each level)	3 x 1 mL
149X	Level Low MiniPak	1 x 3 mL			



**Bio-Rad
Laboratories**

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

**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

Cobas is a trademark of Roche Diagnostics.