



# Liquichek™ D-dimer Control

A Wider Range of Values Across More Platforms



# Liquichek™ D-dimer Control

Provides the laboratory with a liquid, 4 Level, human plasma-based control for monitoring the precision of D-dimer testing on most automated coagulation, chemistry and immunoassay analyzers.



- 3 year shelf life at -20°C to -70°C which allows for fewer crossover studies
- No reconstitution required
- 15 day open-vial stability at 2–8°C

## D-dimer

### Popular Instruments for D-dimer Assays

Instrument	Level Low	Level 1	Level 2	Level 3
Abbott	•	•	•	•
Beckman Coulter® Olympus AU		•	•	•
bioMérieux	•	•	•	•
Biosite Triage® Meter Pro		•	•	•
Diagnostica Stago	•	•	•	•
IL ACL	•	•	•	•
Mitsubishi Pathfast®	•	•	•	
Roche/Hitachi® Cobas	•	•	•	•
Siemens BCS®	•	•	•	•
Siemens Immulite® 2000		•	•	•
Siemens Stratus®	•	•	•	•
Sysmex® CA Series	•	•	•	•

### Target Level

	D-dimer ng/mL (FEU)*
Level Low	250
Level 1	550
Level 2	1600
Level 3	3200

The values listed are approximate targets. The actual observed values will vary depending upon lot number, analyzer, reagent, method, and calibrator.

\* Fibrinogen Equivalent Units (FEU) = twice the mass of the D-DU (D-dimer units).

These are recommended levels. Refer to package insert for complete instrument listing.

## Ordering Information

Cat #	Liquichek™ D-dimer Control	Quantity
27100	Level Low	6 x 1 mL
27101	Level 1	6 x 1 mL
27102	Level 2	6 x 1 mL
27103	Level 3	6 x 1 mL
27102X	Four Level MiniPak (1 of each level)	4 x 1 mL



This control is featured in the Unity™ Interlaboratory Program. Participation in the Unity™ Interlaboratory Program can help improve the quality of your laboratory test results 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 the Bio-Rad office nearest you or visit our website at [www.bio-rad.com/coagcontrols](http://www.bio-rad.com/coagcontrols)*

**Clinical Diagnostics Group**

**Website** [www.bio-rad.com/qualitycontrol](http://www.bio-rad.com/qualitycontrol) **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** 86-21-61698500 **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** 026-674-55-05/06 **Taiwan** 886-2-2578-7189 **Thailand** 662-651-8311 **United Kingdom** +44-(0)20-8328-2000