



Lyphochek[®] Coagulation Control



Lyphochek® Coagulation Control

A human plasma control for monitoring the precision of routine coagulation tests. It has an extended reconstituted stability that helps reduce waste and save money.

- 3 year shelf life at 2–8°C
- 48 hour open-vial stability at 2–25°C
- Representative PT values in seconds and International Normalized Ratio (INR)
- Normal and abnormal levels of PT, APTT, Fibrinogen, Antithrombin III and Thrombin Time
- For use on a variety of systems, including Siemens BCS®, Sysmex® CA, Diagnostica Stago STA®, Instrumentation Laboratory ACL® and ELECTRA™, and Trinity Biotech AMAX



Analytes

Activated Partial Thromboplastin Time (APTT)
Antithrombin III (AT III)
Fibrinogen

Prothrombin Time (PT)
Thrombin Time (TT)

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

Ordering Information

Cat #	Lyphochek® Coagulation Control	Quantity
744	Level 1	12 x 1 mL
745	Level 2	12 x 1 mL
746	Level 3	12 x 1 mL
745X	Trilevel MiniPak (1 of each level)	3 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.



**Bio-Rad
Laboratories**

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

**Clinical
Diagnostics Group**

Website 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** +36-1-459-6100 **India** 1-800-180-1224 **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** 41-61-717-95-55 **Thailand** 662-651-8311 **United Kingdom** +44-(0)20-8328-2000