



Lyphocheck® Endocrine Control



Lyphochek® Endocrine Control

A human serum based control designed to monitor the precision of plasma catecholamine testing procedures by HPLC. The analytes are spiked at critical concentrations to cover the dynamic range of the assays.

- Human serum based
- 2 year shelf life at 2–8°C means fewer crossovers and long term QC monitoring
- 3 day open-vial stability at 2–8°C reduces waste and saves money



Analytes

Dopamine
Epinephrine
Norepinephrine

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

Ordering Information

Cat #	Lyphochek® Endocrine Control	Quantity
535	Bilevel (6 of each level)	12 x 10 mL
535X	Bilevel MiniPak (1 of each level)	2 x 10 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/specialtycontrols

*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** 91-124-4029300 **Israel** 972-3-9636050
Italy +39-02-216091 **Japan** 81-3-6361-7070 **Korea** 82-2-3473-4460 **Mexico** 52(55)5200-0520 **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