



Lyphochek® Urine Bone Markers Control



Lyphochek® Urine Bone Markers Control

Lyphochek® Urine Bone Markers Control is a bilevel, human urine based, third party control designed to monitor the test precision for urinary markers of osteoporosis management and bone resorption.



- Human urine based
- 3 year shelf life at 2–8°C
- 14 day open-vial stability at 2–8°C

Analytes

β_2 Microglobulin*	Deoxypyridinoline (DPD)	Pyridinoline (PYD)*
Creatinine	N-Telopeptide (NTx)	
C-Telopeptide (CTx)	Pyridinium Crosslinks (DPD + PYD)	

* No claims are made for expected values or stability
Refer to the package insert of currently available lots for specific analyte and stability claims

Ordering Information

Cat #	Lyphochek® Urine Bone Markers Control	Quantity
406	Bilevel (3 of each level).....	6 x 2 mL
406X	Bilevel MiniPak (1 of each level).....	2 x 2 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

**Bio-Rad
Laboratories**

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

*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