



# Lyphocheck® Controls for Trace Metals Testing

**BIO-RAD**

## Lyphochek® Urine Metals Control

Incorporates two levels of critical trace elements, heavy metals and organic metabolites, and is ideal for environmental and industrial testing.

- Human urine based
- 3 year shelf life at 2–8°C
- 5 day reconstituted stability at 2–8°C
- Contains 23 trace elements, heavy metals and organic metabolites
- Use with Atomic Absorption, GC-MS, HPLC and ICP

### Analytes

5-Aminolevulinic Acid	Mandelic Acid
Aluminum	Manganese
Antimony	Mercury
Arsenic	Nickel
Cadmium	Pentachlorophenol
Chromium	Phenol
Cobalt	Phenylglyoxylic Acid
Copper	Selenium
Fluoride	Thallium
Hippuric Acid	Trichloroacetic Acid
Lead	Zinc
Magnesium	

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

### Ordering Information

Cat #	Lyphochek® Urine Metals Control	Quantity
400	Level 1 . . . . .	10 x 25 mL
405	Level 2 . . . . .	10 x 25 mL

## Lyphochek® Whole Blood Metals Control

Contains the major heavy metals of toxicological and environmental importance for whole blood analysis.

- Human whole blood based
- 3 year shelf life at 2–8°C
- 14 day reconstituted stability at 2–8°C
- Lead assay values for Atomic Absorption, Anodic Stripping Voltammetry (ASV), Inductively Coupled Plasma-Mass Spectrometry (ICP-MS) and Isotope Dilution Mass Spectrometry (IDMS)

### Analytes

Arsenic
Cadmium
Lead
Mercury
Thallium

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

### Ordering Information

Cat #	Lyphochek® Whole Blood Metals Control	Quantity
527	Level 1 . . . . .	6 x 2 mL
528	Level 2 . . . . .	6 x 2 mL
529	Level 3 . . . . .	6 x 2 mL
528X	Trilevel MiniPak (1 of each level) . . . . .	3 x 2 mL



These controls are 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**

**Bio-Rad  
Laboratories**

**For further information, please contact the Bio-Rad office nearest you or visit our website at [www.bio-rad.com/qualitycontrol](http://www.bio-rad.com/qualitycontrol)**

*Clinical  
Diagnostics Group*

**Website** [www.bio-rad.com/qualitycontrol](http://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