



# Specialty Drug Testing Quality Controls





## QC Solutions for the Unique Challenges of Specialty Drug Testing

**Bio-Rad offers over 300 quality controls, including a comprehensive line of products to monitor drug testing precision. QC solutions are available for the unique challenges of specialty drug testing, such as finding a drug-free baseline serum or a control suitable for the special needs of HPLC testing.**

Along with the right controls for your specialty drug testing program, participation in the Unity™ Interlaboratory Program allows you to compare your results with peer group data from other laboratories. Designed to fit your specific needs, a choice of QC data management solutions offers desktop or web-based alternatives, with QC statistical rules ranging from basic to advanced.



### Lyphochek® Benzo/TCA Control-Set A

Ideal for monitoring the precision of HPLC methods for identification and quantitation of benzodiazepines and tricyclic antidepressant drugs.

- Human serum based
- 3 year shelf life at 2–8°C
- 30 day reconstituted stability at 2–8°C

#### Analytes

Amitriptyline	Diazepam	Nordoxepin
Chlordiazepoxide	Doxepin	Nortriptyline
Desalkylflurazepam	Imipramine	Oxazepam
Desipramine	Nordiazepam	

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



### Lyphochek® Benzo/TCA Control-Set B

Ideal for monitoring the precision of HPLC methods for identification and quantitation of benzodiazepines and antidepressant drugs.

- Human serum based
- 3 year shelf life at 2–8°C
- 14 day reconstituted stability at 2–8°C

#### Analytes

Alprazolam	Norclomipramine
Clomipramine	Trazodone

Clonazepam has been added to this product as a service to laboratories that wish to investigate this analyte.

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



### Lyphochek® Drug Free Serum

Lyophilized serum tested negative for more than 60 therapeutic drugs and illicit compounds.\*

- Drug-free baseline reference for HPLC analysis
- For use as a negative control
- 3 year shelf life at 2–8°C
- 10 day reconstituted stability at 2–8°C

#### Analytes

Acetaminophen	Desipramine	Meperidine	Phenothiazines
Alprazolam	Desmethyldoxepin	Methadone	Phenytol
Amikacin	Diazepam	Methamphetamine	Primidone
Amitriptyline	Digitoxin	Methaqualone	Procainamide
Amphetamine	Digoxin	Methotrexate	Propoxyphene
Caffeine	Disopyramide	Methyprylon	Propranolol
Cannabinoids	Doxepin	Morphine	Pyrimidine
Carbamazepine	Ethosuximide	N-Acetylprocainamide	Quinidine
Chloramphenicol	Flecainide	Netilmicin	Quinine
Chlordiazepoxide	Gentamicin	Nicotine	Salicylate
Cimetidine	Glutethimide	Norclomipramine	Secobarbital
Clomipramine	Haloperidol	Nordiazepam	Theophylline
Clonazepam	Hydromorphone	Nordoxepin	Tobramycin
Cocaine	Ibuprofen	Nortriptyline	Trazodone
Codeine	Imipramine	Oxazepam	Valproic Acid
Cyclosporine	Lidocaine	Phencyclidine	Vancomycin
Desalkylflurazepam	Lithium	Phenobarbital	

\*Refer to the package insert of currently available lots for specific analyte and stability claims, and detection limits of methods used.

## Ordering Information

Cat #	Product Name	Quantity
<b>Lyphochek® Benzo/TCA Control–Set A</b>		
661	Level 1	6 x 5 mL
662	Level 2	6 x 5 mL
660X	Bilevel MiniPak (1 of each level)	2 x 5 mL

Cat #	Product Name	Quantity
<b>Lyphochek® Benzo/TCA Control–Set B</b>		
663	Level 1	6 x 5 mL
664	Level 2	6 x 5 mL
665X	Bilevel MiniPak (1 of each level)	2 x 5 mL

Cat #	Product Name	Quantity
<b>Lyphochek® Drug Free Serum</b>		
456	Single Level	12 x 10 mL
456X	MiniPak	1 x 10 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  
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** 36-1-455-8800 **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)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** 65-6415-3188 **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