



# Lyphocheck® Therapeutic Drug Monitoring Control (TDM)



## Lyphochek® Therapeutic Drug Monitoring Control (TDM)

For labs that run a high volume of serum therapeutic drug tests or prefer dedicated TDM controls. Contains over 40 analytes including markers for monitoring thyroid function.

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



### Analytes

Acetaminophen	Disopyramide	Nortriptyline	T4 (Total)
Amikacin	Estriol (Total)*	Phenobarbital	Theophylline
Amiodarone*	Ethosuximide	Phenytoin	Tobramycin
Amitriptyline	Flecainide*	Phenytoin (Free)	Tricyclic Antidepressants (TCA) Screen
Caffeine	Gentamicin	Primidone	TSH
Carbamazepine	Haloperidol*	Procainamide	Valproic Acid
Carbamazepine (Free)	Imipramine	Propranolol	Valproic Acid (Free)
Chloramphenicol	Lidocaine	Quinidine	Vancomycin
Cortisol	Lithium	Salicylate	
Cyclosporine	Methotrexate	T3 (Total)	
Desipramine	N-Acetylprocainamide (NAPA)	T3 Uptake/T-Uptake	
Digoxin	Netilmicin*	T4 (Free)	

\*No claims are made regarding performance or stability. Refer to myinserts.com or the package insert of currently available lots for specific analyte and stability claims.

### Ordering Information

Cat #	Lyphochek® Therapeutic Drug Monitoring Control (TDM)	Quantity
450	Trilevel (4 of each level) . . . . .	12 x 5 mL
451	Level 1 . . . . .	12 x 5 mL
452	Level 2 . . . . .	12 x 5 mL
453	Level 3 . . . . .	12 x 5 mL
450X	Trilevel MiniPak (1 of each level) . . . . .	3 x 5 mL

### Bio-Rad is Dedicated to Advancing QC Solutions Worldwide

#### Quality Controls

Broad portfolio of independent controls provides an unbiased assessment of methods and system performance.

#### QC Data Management Solutions

Manage and interpret your QC results using the industry-leading Unity™ Interlaboratory Program and benefit from data generated by users worldwide.

#### Customer Support

Receive the backing of full-service technical and customer support trained to handle your inquiries in a prompt, knowledgeable manner.

#### Education

We offer the educational tools and resources to help improve your ability to produce reliable test results.



**Bio-Rad Laboratories**

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

### Clinical Diagnostics Group

**Website** [www.bio-rad.com/qualitycontrol](http://www.bio-rad.com/qualitycontrol) **Australia** 61-2-9914-2800 **Austria** 43-1-877-8901 **Belgium** +32 (3)710-53-00 **Brazil** +55 (31)3689-6600 **Canada** 1-514-334-4372 **China** 86-21-61698500 **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** 1800-180-1224 **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** +7-495-721-1404 **Singapore** 65-6415-3170 **South Africa** 27-11-442-85-08 **Spain** 34-91-590-5200 **Sweden** +46-8-555-127-00 **Switzerland** +41 (0)26-674-55-05/06 **Taiwan** 886-2-2578-7189 **Thailand** 662-651-8311 **United Kingdom** +44 (0)20-8328-2000