



Lyphocheck[®] Therapeutic Drug Monitoring Control (TDM)

BIO-RAD

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 myeinserts.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

Clinical Diagnostics Group

Website 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