



# Lyphocheck<sup>®</sup> Immunoassay Plus Control



# Lyphochek® Immunoassay Plus Control

Provides the clinical laboratory with the most comprehensive control in the world for monitoring immunoassay and therapeutic drug monitoring procedures, eliminating the need for individual controls. Lyphochek® Immunoassay Plus Control is made from human serum and is suitable for most test methods.



- Human serum based
- Assayed for over 1100 methods
- Comprehensive analyte menu
- Includes therapeutic drugs
- 3 year shelf life at 2–8°C

## Analytes

11-Deoxycortisol	Carbamazepine	Fructosamine	Lidocaine	Somatomedin-C
17- $\alpha$ -Hydroxyprogesterone	Carbamazepine (Free)	FSH	Lithium	T3 (Free)
25-Hydroxy Vitamin D	CEA	Gastrin	N-Acetylprocainamide (NAPA)	T3 (Total)
Acetaminophen	Chloramphenicol	Gentamicin	Netilmicin*	T3 Uptake
ACTH	Cortisol	Glucagon	Nortriptyline	T4 (Free)
Aldosterone	Cyclosporine**	hCG	PAP	T4 (Total)
Alphafetoprotein (AFP)	Desipramine	hCG- $\beta$ Subunit	Phenobarbital	TBG
Amikacin	DHEA	hGH	Phenytoin	Testosterone
Amiodarone*	DHEA Sulfate	Imipramine	Phenytoin (Free)	Testosterone (Free)
Amitriptyline	Digoxin	Immunoglobulin A (IgA)	Primidone	Theophylline
Androstenedione	Disopyramide	Immunoglobulin E (IgE)	Procainamide	Thyroglobulin (Tg)
Angiotensin I	Estradiol	Immunoglobulin G (IgG)	Progesterone	Tobramycin
Anti-Thyroglobulin (Anti-Tg)*	Estriol (Free)	Immunoglobulin M (IgM)	Prolactin	Tricyclic Antidepressants (TCA) Screen
Anti-Thyroperoxidase (Anti-TPO)*	Estriol (Total)*	Immunoreactive	PSA	TSH
C-Peptide	Estrogen (Total)	Trypsinogen (IRT)	PSA (Free)	Valproic Acid
Caffeine	Ethosuximide	Insulin	Quinidine	Valproic Acid (Free)
Calcitonin	Ferritin	Iron	Salicylate	Vancomycin
	Flecainide	Iron (TIBC)	SHBG*	Vitamin B <sub>12</sub>
	Folate	LH		

\*No claims are made regarding performance or stability.

\*\* Values are not provided.

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

## Ordering Information

Cat #	Lyphochek® Immunoassay Plus Control	Qty	Cat #	Lyphochek® Immunoassay Plus Control	Qty
370	Trilevel (4 of each level)	12 x 5 mL	373	Level 3	12 x 5 mL
371	Level 1	12 x 5 mL	370X	Trilevel MiniPak (1 of each level)	3 x 5 mL
372	Level 2	12 x 5 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](http://www.QCNet.com).**



**Bio-Rad  
Laboratories**

*For further information, please contact your local Bio-Rad office or learn more online at [www.bio-rad.com/immunocontrols](http://www.bio-rad.com/immunocontrols)*

**Clinical  
Diagnostics Group**

**Website** [www.bio-rad.com/qualitycontrol](http://www.bio-rad.com/qualitycontrol) **U.S.** 1-800-2BIO-RAD (1-800-224-6723) **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** 1-800-180-1224 **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** 7-495-721-14-04 **Singapore** 65-6415-3170 **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