



Liquichek™ Immunoassay Plus Control





Optimize Your Laboratory's Performance with Bio-Rad Multi-Analyte Immunoassay Controls



Liquechek™ Immunoassay Plus Control

A liquid control designed for convenience in today's busy immunoassay laboratory. With the most comprehensive analyte menu in liquid immunoassay controls, this product can replace two or more separate controls for immunoassay and therapeutic drug monitoring procedures. Liquechek™ Immunoassay Plus Control provides an unbiased assessment of your test system.

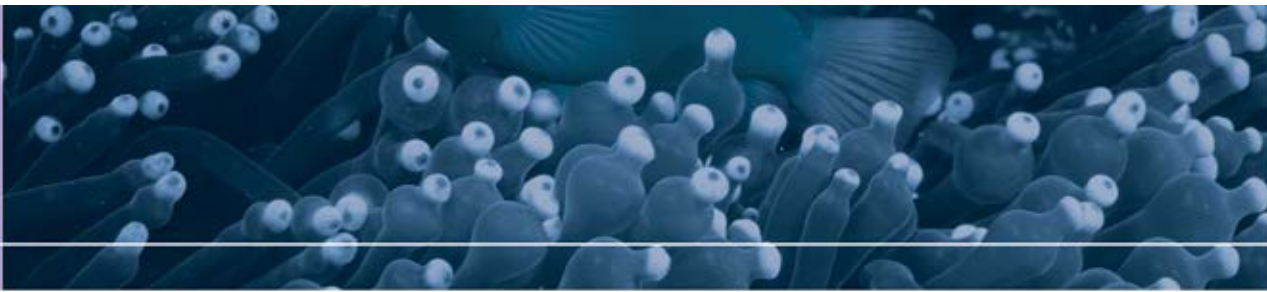
5 mL Fill Size

- Comprehensive IA and TDM control
- Human serum based
- 14 day open-vial stability at 2–8°C for most analytes



2.5 mL Fill Size

- Dimension Vista® vials offer ready-to-use convenience for on-board refrigerated storage
- Offers more than 40 analytes to match the Dimension Vista® immunoassay test menu
- Shared lot numbers and Unity™ peer groups with existing Liquechek™ Immunoassay Plus Control configurations
- 2D barcode technology for quick and easy product insert and vial label scanning
- 2 year shelf life at -20°C to -50°C
- 4 day unopened (unpunctured) stability when stored refrigerated at 2–8°C
- 4 day open-vial (punctured) stability for all analytes when stored at 2–8°C
- Each level sold separately



Analytes for All Fill Sizes

Acetaminophen	DHEA Sulfate ¹	IgM	Salicylate
AFP	Digoxin	Insulin ¹	SHBG ¹
Aldosterone ¹	Disopyramide ¹	Iron	Somatomedin-C* ¹
17-Alpha-Hydroxyprogesterone ¹	Estradiol	Iron (TIBC)* ¹	T3 (Free)
Amikacin	Estriol (Free) ¹	LH	T3 (Total)
Amiodarone* ¹	Estriol (Total)* ¹	Lidocaine	T3 Uptake/T-Uptake
Amitriptyline* ¹	Estrogen (Total) ¹	Lithium	T4 (Free)
Androstenedione ¹	Ethosuximide ¹	NAPA	T4 (Total)
Angiotensin I* ¹	Ferritin	Netilmicin* ¹	TBG ¹
Anti-TG Ab* ¹	Flecainide ¹	Nortriptyline ¹	Testosterone ¹
Anti-TPO Ab* ¹	Folate	PAP ¹	Testosterone (Free) ¹
Caffeine	Fructosamine ¹	Phenobarbital	Theophylline
Carbamazepine	FSH	Phenytoin	Thyroglobulin* ¹
Carbamazepine (Free) ¹	Gentamicin	Phenytoin (Free) ¹	Tobramycin
CEA	hCG ¹	Primidone ¹	Tricyclic Antidepressants (TCA) Screen
Chloramphenicol ¹	hCG-Beta Subunit	Progesterone ¹	TSH
CK-MB Isoenzyme	hGH ¹	Prolactin	Valproic Acid
Cortisol ¹	25-Hydroxy Vitamin D* ¹	Propranolol ¹	Valproic Acid (Free) ¹
Cyclosporine ¹	Ibuprofen ¹	PSA	Vancomycin
11-Deoxycortisol* ¹	Imipramine* ¹	PSA (Free)	Vitamin B ₁₂
Desipramine ¹	IgA	PTH-MM* ¹	
DHEA ¹	IgE	Quinidine ¹	
	IgG		

*No claims are made regarding performance or stability.

¹Not included for use on the Dimension Vista[®] test menu.

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

Ordering Information

Cat #	Product Name	Quantity
Liquichek™ Immunoassay Plus Control		
360	Trilevel (4 of each level)	12 x 5 mL
361	Level 1	12 x 5 mL
362	Level 2	12 x 5 mL
363	Level 3	12 x 5 mL
360X	Trilevel MiniPak (1 of each level)	3 x 5 mL
267	Level 1	12 x 2.5 mL
268	Level 2	12 x 2.5 mL
269	Level 3	12 x 2.5 mL
268X	Trilevel MiniPak (1 of each level)	3 x 2.5 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.



**Bio-Rad
Laboratories**

For further information, please contact the Bio-Rad office nearest you or visit our website at www.bio-rad.com/iacontrols

**Clinical
Diagnostics Group**

Website www.bio-rad.com/qualitycontrol **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** 86-21-61698500 **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** 7-495-721-14-04 **Singapore** 65-6415-3170 **South Africa** 27-11-442-85-08 **Spain** 34-91-590-5200 **Sweden** 46-8-555-127-00 **Switzerland** +41 (0)26-674-52-01/02/03 **Taiwan** 886-2-2578-7189 **Thailand** 662-651-8311 **United Kingdom** +44 (0)20-8328-2000