



# Independent and Complete Control for Immunoassay

Introducing Lyphocheck® Specialty Immunoassay Control with Procalcitonin





## Introducing Lyphochek® Specialty Immunoassay Control with Procalcitonin and Intact PTH

**Complementing an extensive premium menu with even greater specialty coverage**

This new product provides you unbiased independent quality control for a key marker of sepsis. Taken together with intact PTH for monitoring thyroid surgery, this control has utility in Critical Care, Emergency and Operating Room settings.

### Lyphochek® Specialty Immunoassay Control

An assayed control designed to include specialty analytes such as Procalcitonin, Intact PTH, and 25-OH Vitamin D.

- Lyophilized, human serum based
- Three distinct levels
- 3 year shelf life at 2-8°C
- 30 day open-vial stability at 2-8°C for most analytes after reconstitution
- Assayed values available for both automated and manual methods

#### Analytes

25-Hydroxy-Vitamin D  
Erythropoietin (EPO)  
Intact PTH  
Procalcitonin\* **NEW**  
SHBG (Sex Hormone Binding Globulin)

## Or Select Intact PTH and Thyroid Antibodies

**Add just the specialty analytes you need and only when you need them**

### Liquichek™ Specialty Immunoassay Control

An assayed control designed to include specialty analytes such as Anti-Thyroglobulin, Anti-Thyropoxidase, Intact PTH, and 25-OH Vitamin D.

- Liquid, human serum based
- Three distinct levels
- 3 year shelf life at -20°C to -70°C
- 30 day open-vial stability at 2-8°C for most analytes
- Assayed values available for most automated immunoassay analyzers and manual methods
- 25-OH Vitamin D insert listing also includes LC-MS-MS values for Vitamin D2 and D3
- Level LTA designed to provide low levels of Anti-Tg and Anti-TPO values for most platforms.

#### Analytes

25-Hydroxy Vitamin D  
Anti-Thyroglobulin (Anti-Tg)  
Anti-Thyropoxidase (Anti-TPO)  
C-Peptide  
Erythropoietin (EPO)  
Insulin Growth Factor I (IGF-I)  
Intact PTH  
Osteocalcin





## Consolidated Quality Control for Immunoassay and Tumor Markers

**Greater efficiency for your integrated analyzer**

### Liquichek™ Immunoassay Premium Control

Using Liquichek™ Immunoassay Premium Control allows your laboratory to replace two or more individual QC products with a single control while maintaining the same QC program for data review and peer group reports.

By bringing together Immunoassays, Specialty Immunoassays (25-OH Vitamin D, SHBG), Tumor Markers, and Therapeutic Drug Monitoring analytes, Bio-Rad has created an effective way to maximize the power of your quality control plan.

This new premium QC product is an unbiased independent control providing assayed values for most major integrated immunoassay platforms. International Organization for Standardization (ISO) 15189 recommends the use of independent control materials; either instead of, or in addition to, any control materials supplied by the reagent or instrument manufacturers.

- Liquid, human serum based
- Three distinct levels
- 2 year shelf life at -20°C to -70°C
- 14 day open-vial stability at 2-8°C for most analytes
- Includes Tumor Markers

#### Analytes

##### Immunoassay

Acid Phosphatase, Prostatic  
AFP  
Cortisol  
DHEA-Sulfate  
Estradiol  
Estrinol, Free  
Ferritin  
Folate  
FSH  
hCG, Total beta  
hGH  
Insulin  
LH  
Progesterone  
Prolactin  
T3 Uptake/T Uptake  
T3, Free  
T3, Total  
T4, Free  
T4, Total  
TBG  
Testosterone  
TSH  
Vitamin B<sub>12</sub>

##### Immunology

IgA  
IgE  
IgG  
IgM

##### Specialty Immunoassay

25-Hydroxy Vitamin D (Total)  
SHBG (Sex Hormone Binding Globulin)

##### TDM

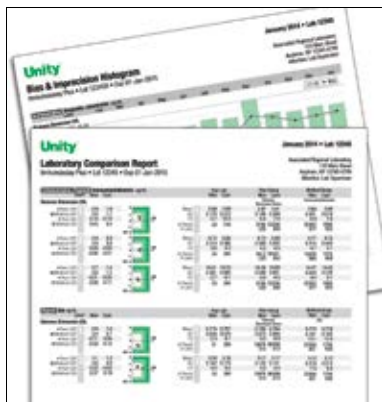
Acetaminophen  
Amikacin  
Caffeine  
Carbamazepine (Tegretol)  
Digoxin  
Disopyramide  
Ethosuximide  
Gentamicin  
Lidocaine  
Lithium  
NAPA (N-Acetyl Procainamide)  
Phenobarbital  
Phenytoin  
Phenytoin, Free  
Primidone  
Procainamide  
Quinidine  
Salicylate  
Theophylline  
Tobramycin  
Tricyclic Antidepressant (TCA) Screen  
Valproic Acid  
Valproic Acid, Free  
Vancomycin

##### Tumor Marker

β-2-Microglobulin  
CA 125  
CA 15-3  
CA 19-9  
CEA (Carcinoembryonic Antigen)  
Cyfra 21-1\*\*  
PSA, Free  
PSA, Total

**Complete your Immunochemistry coverage with Liquid Assayed Multiquel® Premium Control**





## Unity™ QC Data Management Solutions

Bio-Rad Quality Controls are part of the Unity™ Program with its industry-leading features for benchmarking laboratory performance.

- Benefit from peer data generated from more than 40,000 instruments worldwide
- Utilize real-time access to peer group comparisons to help with troubleshooting
- Consolidate your entire laboratory into one QC system
- Facilitate run-time decisions with optimized Westgard rules
- Facilitate compliance with simple tools to review rule violations and document corrective actions
- Define your own quality requirements and track key quality performance indicators

You can consolidate your entire laboratory into one QC system through our Unity™ Program. This sophisticated software includes best-in-class tools to support releasing results, benchmarking laboratory performance, and designing your QC system.

## Ordering Information

Cat #	Product Name	Quantity	Cat #	Product Name	Quantity
<b>Lyphocheck® Specialty Immunoassay Control</b>			<b>Liquichek™ Immunoassay Premium Control</b>		
27124	Level 1 . . . . .	6 x 2 mL	27110	Trilevel (2 of each level) . . . . .	6 x 5 mL
27125	Level 2 . . . . .	6 x 2 mL	27111	Level 1 . . . . .	6 x 5 mL
27126	Level 3 . . . . .	6 x 2 mL	27112	Level 2 . . . . .	6 x 5 mL
27125X	Trilevel MiniPak (1 of each level) . . . . .	3 x 2 mL	27113	Level 3 . . . . .	6 x 5 mL
<b>Liquichek™ Specialty Immunoassay Control</b>			<b>Liquid Assayed Multiquel® Premium Control</b>		
359	Level LTA . . . . .	6 x 5 mL	284	Level 1 . . . . .	6 x 5 mL
364	Level 1 . . . . .	6 x 5 mL	285	Level 2 . . . . .	6 x 5 mL
365	Level 2 . . . . .	6 x 5 mL	286	Level 3 . . . . .	6 x 5 mL
366	Level 3 . . . . .	6 x 5 mL	285X	Trilevel MiniPak (1 of each level) . . . . .	3 x 5 mL
359X	Four Level MiniPak (1 of each level) . . . . .	4 x 5 mL			



**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) **Australia** 61-2-9914-2800 **Austria** 43-1-877-8901 **Belgium** 32-9-385-5511 **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