



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
Estradiol, 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 B₁₂

Immunology

IgA
IgE
IgG
IgM

Specialty Immunoassay

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

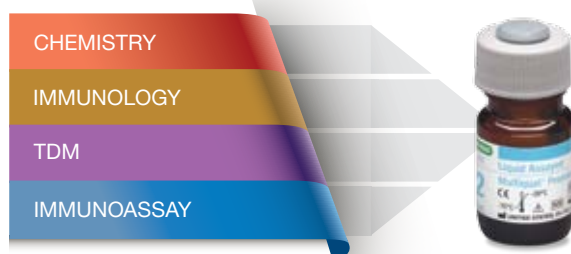
TDM

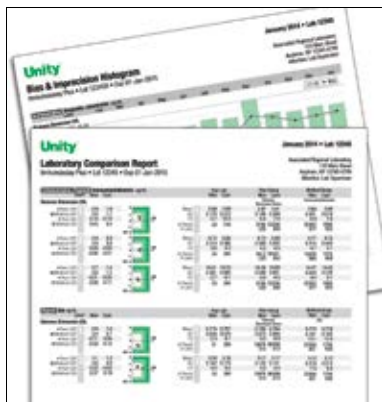
Acetaminophen
Amikacin
Caffeine
Carbamazepine (Tegretol)
Digoxin
Disopyramide
Ethosuximide
Gentamicin
Lidocaine
Lithium
NAPA (N-Acetyl Procainamide)
Phenobarbital
Phenytoin
Phenytoin, Free
Primidone
Procainamide
Quinidine
Salicylate
Theophylline
Tobramycin
Tricyclic Antidepressant (TCA) Screen
Valproic 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
Lyphocheck® Specialty Immunoassay Control			Liquichek™ Immunoassay Premium Control		
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
Liquichek™ Specialty Immunoassay Control			27110X MiniPak (1 of each level)		
359	Level LTA	6 x 5 mL	Liquid Assayed Multiquel® Premium Control		
364	Level 1	6 x 5 mL	284	Level 1	6 x 5 mL
365	Level 2	6 x 5 mL	285	Level 2	6 x 5 mL
366	Level 3	6 x 5 mL	286	Level 3	6 x 5 mL
359X	Four Level MiniPak (1 of each level)	4 x 5 mL	285X	Trilevel MiniPak (1 of each level)	3 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

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** +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