

Intelligent QC Starts with **InteliQ**



Load-and-Go Quality Controls
for Workflow Solutions

BIO-RAD

InteliQ

The Smart QC Choice for Every Lab



Optimize Workflow

Increase productivity and improve turnaround time by eliminating manual steps.



Increase Efficiency

Unique sample tracking and easy XML data upload of lot-specific QC values.



Improve Performance

Access the Unity data management program and large peer groups.



Achieve Quality Goals

Use advanced Unity software tools to support your lab's risk management program and simplify compliance reporting.



Enhance Your Load-and-Go QC with Unity QC Data Management Solutions







InteliQ and Unity

Powerful tools for streamlined workflow efficiency and robust data management

Automate your QC with InteliQ on next generation platforms. These barcoded, load-and-go controls reduce hands-on time and manual errors, streamlining your workflow and giving you more freedom. Together with Unity's advanced data management tools, InteliQ controls improve laboratory efficiency.

InteliQ Quality Controls and Unity software offer you the following benefits:

<p>InteliQ Tubes</p> 	<ul style="list-style-type: none"> ▪ Eliminate manual pour-off with ready-to-use “load and go” convenience to minimize handling errors ▪ Reduce use of consumables and minimize laboratory waste ▪ Conveniently store refrigerated on-board the analyzer*
<p>InteliQ Barcode Labels</p> 	<ul style="list-style-type: none"> ▪ Each InteliQ tube is labeled with a unique sample identifier that provides traceability ▪ Eliminates manual QC order entry and prevents handling errors for a simpler, more efficient workflow
<p>eData</p> 	<ul style="list-style-type: none"> ▪ Import value assignment data to the analyzer with Bio-Rad's proprietary format for embedded data security** ▪ Eliminates manual data input and prevents transcription errors ▪ Provides open-vial stability information for available analytes
<p>Unity</p> 	<ul style="list-style-type: none"> ▪ Access the industry leading Unity data management program and global peer group comparisons for greater confidence in reliability of patient results ▪ Use advanced Unity software tools to support your lab's risk management program and simplify compliance reporting to achieve quality goals

*Optional Atellica Solution.

**Please refer to the package insert for platform-specific stability and performance claims.

InteliQ Quality Controls

InteliQ Assayed Multiquel Control



Comprehensive range of 80 analytes to maximize testing efficiency for automated platforms.

- Liquid, human serum based
- 3 year shelf life at -20°C to -70°C
- 14 day on-board stability at 2-8°C for most analytes*
- 7 day open-vial stability at 2-8°C for most analytes

Analytes

α-1-Antitrypsin	Calcium	Haptoglobin	Salicylate
α-1-Globulin	Calcium (Ionized)	Immunoglobulin A (IgA)	Sodium
α-2-Globulin	Carbamazepine	Immunoglobulin G (IgG)	T3 (Total)
β-Globulin	Carbon Dioxide (CO ₂)	Immunoglobulin M (IgM)	T3 Uptake/T Uptake
γ-Globulin	Ceruloplasmin	Iron	T4 (Free)
γ-Glutamyltransferase (GGT)	Chloride	Iron (TIBC)	T4 (Total)
Acetaminophen	Cholesterol (HDL)	Iron (UIBC)	Theophylline
Acid Phosphatase (Total)	Cholesterol (LDL)	Lactate (Lactic Acid)	Thyroid Stimulating Hormone (TSH)
Albumin	Cholesterol (Total)	Lactate Dehydrogenase (LDH)	Tobramycin
Albumin-to-Globulin Ratio	Cholinesterase	LAP Arylamidase	Transferrin
Alkaline Phosphatase	Complement C3	Lipase	Triglycerides
ALT/SGPT	Complement C4	Lithium	Urea
Amikacin	Copper	Magnesium	Urea Nitrogen (BUN)
Amylase	Cortisol	Osmolality	Uric Acid
Amylase (Pancreatic)	Creatine Kinase (CK)	Phenobarbital	Valproic Acid
Apolipoprotein A-I	Creatinine	Phenytoin	Vitamin B ₁₂
Apolipoprotein B	Digoxin	Phospholipids	Zinc
AST/SGOT	Ethanol	Phosphorus	
Bilirubin (Direct)	Ferritin	Potassium	
Bilirubin (Neonatal)	Gentamicin	Prealbumin	
Bilirubin (Total)	Glucose	Protein Serum (Total)	

InteliQ Cardiac Markers Plus Control LT



Multi-analyte control designed to monitor precision of cardiac assessment assays.

- Liquid, human serum based
- 3 year shelf life at -20°C to -70°C
- 20 day on-board stability at 2-8°C for most analytes*
- 10 day open-vial stability at 2-8°C for most analytes

NEW Analyte

Analytes

BNP	Myoglobin
CK (Total)	NT-proBNP
CK-MB Isoenzyme	Troponin I**
CRP	Troponin T
Digitoxin	

InteliQ Diabetes Control



Trilevel HbA1c control designed to monitor the precision of test procedures associated with diabetes monitoring for Hemoglobin A1C.

- Liquid, human whole blood based
- 3 year shelf life at -20°C to -70°C
- 45 day on-board stability at 2-8°C*
- 10 day open-vial stability at 2-8°C

Analytes

Hemoglobin A1C

*Please refer to the package insert for platform-specific stability and performance claims.

**High-sensitivity Troponin I available.

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



InteliQ Ethanol/Ammonia Control



Normal, elevated and toxic concentrations of Ethanol and Ammonia to monitor precision across the clinical range.

- 2 year shelf life at 2-8°C
- Ammonia:
 - 13 day on-board stability at 2-8°C*
 - 4 day open-vial stability at 2-8°C
- Ethanol:
 - 3 day on-board stability at 2-8°C*
 - 3 day open-vial stability at 2-8°C

Analytes

Ammonia
Ethanol

InteliQ Homocysteine Control



Monitors precision of Homocysteine assays.

- Liquid, human serum based
- 3 year shelf life at -20°C to -70°C
- 30 day on-board stability at 2-8°C*
- 5 day open-vial stability at 2-8°C

Analyte

Homocysteine

InteliQ Immunoassay Plus Control



Offers a wide range of popular immunoassay analytes. Provides an efficient solution for laboratories that focus on routine testing.

- Liquid human serum based
- 28 month shelf life at -20°C to -70°C
- 14 day on-board stability at 2-8°C*
- 10 day open-vial stability at 2-8°C

NEW Analytes

Analytes

Acetaminophen
AFP
Amikacin
Androstenedione
Caffeine
Carbamazepine
CEA
CK-MB Isoenzyme
Cortisol
Dehydroepiandrosterone-Sulfate
(DHEA-Sulfate)
Digoxin
Estradiol
Estriol (Free)

Ferritin
Folate
FSH
Gentamicin
hCG
hGH
Immunoglobulin A (IgA)
Immunoglobulin E (IgE)
Immunoglobulin G (IgG)
Immunoglobulin M (IgM)
Insulin
Iron
Lidocaine
Lithium

Luteinizing Hormone (LH)
N-Acetyl Procainamide (NAPA)
Phenobarbital
Phenytoin
Primidone
Procainamide
Progesterone
Prolactin
PSA (Free)
PSA (Total)
Quinidine
Salicylate
SHBG
T3 (Free)

T3 (Total)
T3 Uptake/T Uptake
T4 (Free)
T4 (Total)
Testosterone
Theophylline
Thyroid Stimulating Hormone (TSH)
Thyroxine Binding Globulin (TBG)
Tobramycin
Tricyclic Antidepressants (TCA)
Valproic Acid
Vancomycin
Vitamin B₁₂

*Please refer to the package insert for platform-specific stability and performance claims. Refer to myelinerts.com or the package insert of currently available lots for specific analyte and stability claims.



IntelIQ Immunology Control



Comprehensive menu of serum proteins and analytes, including Rheumatoid Factor.

- Liquid human serum based
- 2 year shelf life at -20°C to -70°C
- 30 day on-board stability*
- 10 day open-vial stability at 2-8°C

NEW Analytes

Analytes

α -1-Acid Glycoprotein (AAG)
 α -1-Antitrypsin (AAT)
 α -2-Macroglobulin
 β -2-Microglobulin
 ADNase B
 Albumin
 Antistreptolysin O (ASO)
 Anti-Cyclic Citrullinated Peptide
 (Anti-CCP)

Antithrombin III
 Apolipoprotein A-1
 Apolipoprotein B
 C1 Inhibitor
 Ceruloplasmin
 Complement C3
 Complement C4
 C-Reactive Protein (CRP)
 Cystatin C

Ferritin
 Haptoglobin
 Immunoglobulin A (IgA)
 Immunoglobulin E (IgE)
 Immunoglobulin G (IgG)
 Immunoglobulin G Subclasses 1-4
 Immunoglobulin M (IgM)
 Kappa Light Chain
 Lambda Light Chain

Lipoprotein (a)
 Prealbumin
 Properdin Factor B
 Protein Serum (Total)
 Retinol Binding Protein (RBP)
 Rheumatoid Factor
 Transferrin

IntelIQ Specialty Immunoassay Control



Multi-analyte control for monitoring the precision of specialty immunoassay testing.

- Liquid, human serum based
- 2 year shelf life at -20°C to -70°C
- 30 day on-board stability at 2-8°C*
- 10 day open-vial stability at 2-8°C

Analytes

25-Hydroxy Vitamin D
 Anti-Thyroglobulin (Anti-Tg)
 Anti-Thyropoxidase (Anti-TPO)
 C-Peptide

Erythropoietin (EPO)
 Insulin Growth Factor I (IGF-I)/Somatomedin C
 Intact PTH
 Osteocalcin

InteliQ Tumor Marker Control



Comprehensive, assayed, multi-analyte control for monitoring precision of immunoassay-based oncology testing procedures.

- Liquid, human serum based
- 28 month shelf life at -20°C to -70°C
- 30 day on-board stability at 2-8°C*
- 10 day open-vial stability at 2-8°C

Analytes

α-Fetoprotein (AFP)
β-2-Microglobulin
CA 15-3
CA 19-9
CA 27.29
CA 72-4**
CA 125
Carcinoembryonic Antigen (CEA)
Cyfra 21-1**
Ferritin
hCG

HE4
Her-2/neu
Insulin Growth Factor I (IGF-I)
PAP
ProGRP**
Prolactin
PSA (Free)
PSA (Total)
SCC**
Thyroglobulin (Tg)

**International Use Only

InteliQ Urine Chemistry Control



Liquid, bilevel human urine based control ideal for monitoring urine chemistry test procedures.

- Offers 19 analytes
- 2 year shelf life at 2-8°C
- 30 day on-board stability at 2-8°C*
- 10 day open-vial stability at 2-8°C

Analytes

Amylase
Calcium
Chloride
Cortisol
Creatinine

Glucose
Magnesium
Microalbumin
Osmolality
pH

Phosphorus
Potassium
Pregnancy/hCG (Qualitative)
Protein (Total)
Sodium

Specific Gravity
Urea
Urea Nitrogen (BUN)
Uric Acid



XML value assignment available for next generation platforms such as the Siemens Atellica Solution.

Easily upload product parameters and value assignment data to your Atellica Solution using eInserts XML, a standard data exchange format that saves manual data input time, while reducing the risk of transcription errors.

Please refer to the eInserts XML flyer for step-by-step instructions.

Visit www.myinserts.com for easy access to your eData



Ordering Information

Cat # Description

InteliQ Assayed Multiquel Control

12008256	Level 1.	12 x 3 mL
12008257	Level 2.	12 x 3 mL
12008258	Level 3.	12 x 3 mL
12008259	Trilevel MiniPak (1 of each level).	3 x 3 mL

InteliQ Cardiac Markers Plus Control LT

12009955	Level 1.	6 x 3 mL
12009958	Level 1A.	6 x 3 mL
12009959	Level 1B	6 x 3 mL
12009960	Level 1C	6 x 3 mL
12009956	Level 2.	6 x 3 mL
12009957	Level 3.	6 x 3 mL
12009961	Level 1 MiniPak	1 x 3 mL
12009964	Level 1A MiniPak	1 x 3 mL
12009965	Level 1B MiniPak	1 x 3 mL
12009966	Level 1C MiniPak	1 x 3 mL
12009962	Level 2 MiniPak	1 x 3 mL
12009963	Level 3 MiniPak	1 x 3 mL

InteliQ Diabetes Control

12008309	Level 1.	6 x 2 mL
12008310	Level 2.	6 x 2 mL
12008311	Level 3.	6 x 2 mL
12008312	Trilevel MiniPak (1 of each level).	3 x 2 mL

InteliQ Ethanol/Ammonia Control

12008299	Level 1.	6 x 2 mL
12008300	Level 2.	6 x 2 mL
12008301	Level 3.	6 x 2 mL
12008302	Trilevel MiniPak (1 of each level).	3 x 2 mL

InteliQ Homocysteine Control

12004309	Level 1.	6 x 1 mL
12004310	Level 2.	6 x 1 mL
12004311	Level 3.	6 x 1 mL
12004312	Trilevel MiniPak (1 of each level).	3 x 1 mL

Cat # Description

InteliQ Immunoassay Plus Control

12009948	Level 1.	12 x 4 mL
12009949	Level 2.	12 x 4 mL
12009950	Level 3.	12 x 4 mL
12009951	Trilevel MiniPak (1 of each level).	3 x 4 mL

InteliQ Immunology Control

12009941	Level 1.	6 x 3 mL
12009942	Level 2.	6 x 3 mL
12009943	Level 3.	6 x 3 mL
12009944	Trilevel MiniPak (1 of each level).	3 x 3 mL

InteliQ Specialty Immunoassay Control

12008276	Level LTA.	6 x 4 mL
12008277	Level 1.	6 x 4 mL
12008278	Level 2.	6 x 4 mL
12008279	Level 3.	6 x 4 mL
12008280	Four Level MiniPak (1 of each level).	4 x 4 mL

InteliQ Tumor Marker Control

12008289	Level 1.	6 x 2 mL
12008290	Level 2.	6 x 2 mL
12008291	Level 3.	6 x 2 mL
12008292	Trilevel MiniPak (1 of each level).	3 x 2 mL

InteliQ Urine Chemistry Control

12009995	Level 1.	12 x 3 mL
12009996	Level 2.	12 x 3 mL
12009997	Bilevel MiniPak (1 of each level).	2 x 3 mL

Bio-Rad packaging is certified environmentally friendly.

EQAS

An independent, external assessment of performance in comparison to your peers.

QCNet.com/eqas

Independent QC

Ongoing, proactive, unbiased daily QC that helps identify errors as they occur or begin to trend.

QCNet.com/independentqc

Unity

QC Data Management tools that help you create a strategy to reduce risk and streamline QC workflow.

QCNet.com/datamanagement



**Bio-Rad
Laboratories, Inc.**

For further information, please contact the Bio-Rad office nearest you or visit our website at <https://qcnet.com/inteliq/index.html>

Clinical
Diagnostics Group

Website www.bio-rad.com/diagnostics Australia +61 (2) 9914 2800 Austria +43 (0) 1 877 89 01 9 Belgium +32 (0) 3 710 53 00 Brazil +55 11 3065 7550 Canada +1 514 334 4372
 China +86 21 6169 8500 Czech Republic +420 241 431 660 Denmark +45 44 52 10 00 Finland +358 9 804 22 00 France +33 (0)1 47 95 60 00 Germany +49 (0) 89 31884 393
 Greece +30 210 7774396 Hong Kong +85 2 2789 3300 Hungary +36 1 459 6190 India +91 124 4029300 Israel +972 03 963 6025 Italy +39 024 94 86 600
 Japan +81 3 6361 7070 Korea +82 080 007 7373 Mexico +52 (55) 5488 7670 The Netherlands +31 (0)318 540 666 New Zealand +64 (9)415 2280 Norway +47 23 38 41 30
 Poland +48 22 331 99 99 Portugal +351 21 47 27 700 Russia +7 495 721 1404 Singapore +65 6415 3170 South Africa +27 11 442 8508 Spain +34 91 490 6580
 Sweden +46 844 98053 Switzerland +41 (0) 61 717 9555 Taiwan +886 (2) 2578-7189 Thailand (662) 651 8311 United Kingdom +44 (0)1923 471301