

Quality Controls for Workflow Solutions

InteliQ Analyte Chart



- Analyte available
- Analyte present in endogenous levels

Test Menu	Analytes	InteliQ Assayed Multiquant	InteliQ Cardiac Markers Plus LT	InteliQ Diabetes	InteliQ Ethanol/Ammonia	InteliQ Homocysteine	InteliQ Immunoassay Plus	InteliQ Immunology	InteliQ Specialty Immunoassay	InteliQ Tumor Marker	InteliQ Urine Chemistry
Anemia	Ceruloplasmin	○						●			
	Erythropoietin (EPO)								●		
	Ferritin	○					●	●		●	
	Folate						●				
	Vitamin B ₁₂	○					●				
Bone Metabolism	25-Hydroxy Vitamin D								●		
	Intact PTH								●		
	Osteocalcin								●		
Cancer/Oncology	α-Fetoprotein (AFP)						●			●	
	CA 15-3									●	
	CA 19-9									●	
	CA 27.29									●	
	CA 72-4*									●	
	CA 125									●	
	Carcinoembryonic Antigen (CEA)						●			●	
	Cyfra 21-1*									●	
	HE4									●	
	HER-2/neu									●	
	PAP									●	
	ProGRP*									●	
	PSA (Free)							●		●	
	PSA (Total)							●		●	
SCC*									●		
Thyroglobulin									●		
Cardiac	BNP		●								
	CK (Total)		●								
	CK-MB Isoenzyme		●				●				
	CRP		●								
	Myoglobin		●								
	NT-proBNP		●								
	Troponin I**		●								
Troponin T		●									
Diabetes	C-Peptide								●		
	Hemoglobin A1C			●							
	Insulin						●				
Fertility/Pregnancy	Androstenedione						●				
	DHEA-S						●				
	Estradiol						●				
	Estriol (Free)						●				
	FSH						●				
	hCG (Total)						●			●	
	LH						●				
	Pregnancy/hCG (Qualitative)										●
	Progesterone						●				
	Prolactin						●			●	
SHBG (Sex Hormone Binding Globulin)						○					
Testosterone						●					

* International use only.

** Available high-sensitivity Troponin I.



Quality Controls for Workflow Solutions

InteliQ Analyte Chart



- Analyte available
- Analyte present in endogenous levels

Test Menu

Analytes

		InteliQ Assayed Multiquant	InteliQ Cardiac Markers Plus LT	InteliQ Diabetes	InteliQ Ethanol/Ammonia	InteliQ Homocysteine	InteliQ Immunoassay Plus	InteliQ Immunology	InteliQ Speciality Immunoassay	InteliQ Tumor Marker	InteliQ Urine Chemistry
Specific Proteins	α-1-Acid Glycoprotein (AAG)							●			
	α-1-Antitrypsin (AAT)	○						●			
	α-2-Macroglobulin							●			
	ADNase B							●			
	Anti-Cyclic Citrullinated Peptide (Anti-CCP)							●			
	Antithrombin III (AT III)							●			
	Apolipoprotein A-1	○						●			
	Apolipoprotein B	○						●			
	β-2-Microglobulin							●		●	
	C1 Inhibitor							●			
	Complement C3	○						●			
	Complement C4	○						●			
	Cystatin C							○			
	Haptoglobin	○						●			
	IgG Subclass 1							●			
	IgG Subclass 2							●			
	IgG Subclass 3							●			
	IgG Subclass 4							●			
	Immunoglobulin A (IgA)	○						○	●		
	Immunoglobulin G (IgG)	○						○	●		
	Immunoglobulin M (IgM)	○						○	●		
	Kappa Light Chain							●			
	Lambda Light Chain							●			
	Lipoprotein (a)							●			
	pH										●
	Prealbumin	○						●			
	Propterin Factor B							●			
	Protein Serum (Total)	●						●			
	Retinol Binding Protein (RBP)							●			
	Rheumatoid Factor							●			
	Specific Gravity										●
	Transferrin	○						●			



Bio-Rad
Laboratories, Inc.

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

Clinical
Diagnostics Group

Website www.bio-rad.com/qualitycontrol **U.S.** 1 800 224 6723 **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

BIO-RAD and INTELIQ are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. Atellica is a registered trademark of Siemens Healthineers.

© 2021 Bio-Rad Laboratories, Inc. Printed in USA QSD21-0102 Q-1228 Rev. 04/2021

