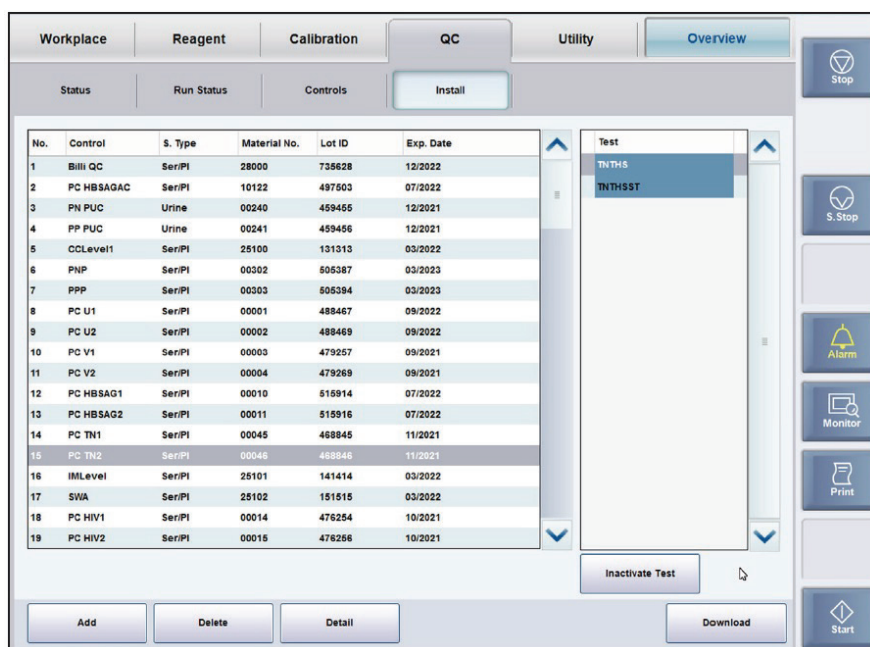


# InteliQ Quick Reference Guide for Roche Cobas 8000 System (w/o API functionality)

## 1. QC Set-Up (Add controls manually)

1. Select a blank line from the control list on the **QC>Install** screen.
2. Choose **Add**.



3. Type all necessary data in the **ADD Control** window.

The 'Add Control' dialog box is shown with the following fields filled out:

- Control: Non-Roche
- Sample Type: Ser/PI
- Material No.: 001
- Lot No.: 123456
- Expiration: 12 - 12

Buttons for 'Cancel' and 'OK' are visible at the bottom.

4. Choose **OK**. The new control appears on the control list

To run control tests with the newly added controls, it is necessary to manually enter the target values for each test and then to activate these tests.

5. Choose **Edit** to enter target mean, target SD, and confidence limits if necessary.

Test	Reagent Lot No.	Individual		Cumulative		Confidence	
		T.Mean	T.SD	T.Mean	T.SD	Low	High
ALB2		47.6	2.9	47.6	2.9	44.7	50.5
ALTL		50.1	3.0	50.1	3.0	47.1	53.1
ASTL		47.6	2.9	47.6	2.9	44.7	50.5
ASTLP		48.9	2.9	48.9	2.9	46.0	51.8
BILTS		1.14	0.07	1.14	0.07	1.07	1.21
CA		2.09	0.08	2.09	0.08	2.01	2.17
CREJ2		96.1	5.8	96.1	5.8	90.3	101.9
CI		83.9	2.5	83.9	2.5	78.9	88.9
GLUC2		5.10	0.25	5.10	0.25	4.85	5.35
GLUC3		5.10	0.25	5.10	0.25	4.85	5.35
IGG-2		7.07	0.42	7.07	0.42	6.65	7.49
K		3.41	0.10	3.41	0.10	3.21	3.61
Na		121	4	121	4	113	129
SBIL5		19.5	1.2	19.5	1.2	18.3	20.7

6. Choose **Active Test** to activate the tests for the new control.

- On the **QC>Install**, choose the appropriate control name in the list box on the left side of the screen.
- Choose all tests to be activated in the list on the right side of the screen.
- Choose **Activate Test**. The test name is highlighted when the test is activated.

Status	Run Status	Individual	Cumulative	Control	Install			
Control	S. Type	QC Number	Lot No.	Test	Individual T.Mean T.SD	Cumulative T.Mean T.SD	Confidence Low High	
1	PNP	Ser PI	302	15654700	ALB2	47.6 2.9	47.6 2.9	44.7 50.5
2	PFU	Ser PI	301	15041500	ALTL	50.1 3.0	50.1 3.0	47.1 53.1
3	PNU	Ser PI	300	15040300	ASTL	47.6 2.9	47.6 2.9	44.7 50.5
4	PPP	Ser PI	303	15654800	ASTLP	48.9 2.9	48.9 2.9	46.0 51.8
5	PCAIN	Suprnt	208	62936900	BILTS	1.14 0.07	1.14 0.07	1.07 1.21
6	PBIL	Ser PI	306	15724500	CA	2.09 0.08	2.09 0.08	2.01 2.17
7	HBCN	Suprnt	357	62341400	CREJ2	96.1 5.8	96.1 5.8	90.3 101.9
8	HBCP	Suprnt	358	61849900	CI	83.9 2.5	83.9 2.5	78.9 88.9
9	PC CARD1	Ser PI	47	16011600	GLUC2	5.10 0.25	5.10 0.25	4.85 5.35
10	PC CARD2	Ser PI	48	16011700	GLUC3	5.10 0.25	5.10 0.25	4.85 5.35
11	PC THYRO1	Ser PI	52	15834200	IGG-2	7.07 0.42	7.07 0.42	6.65 7.49
12	PC THYRO2	Ser PI	53	15804700	K	3.41 0.10	3.41 0.10	3.21 3.61
13	PC A1	Ser PI	54	15390900	Na	121 4	121 4	113 129
14	PC A2	Ser PI	55	15221300	SBIL5	19.5 1.2	19.5 1.2	18.3 20.7
15	PC A3	Ser PI	56	15221400	TP2	64.0 2.6	64.0 2.6	61.4 66.6
16	PC TM1	Ser PI	6	15428200	TRSP2	1.95 0.12	1.95 0.12	1.83 2.07
17	PC TM2	Ser PI	7	15428300	wtzecal			
18	biorad	Ser PI	802	0101010				

## 7. Assign a control to Rack

1. On **QC>Control** choose **Rack Assignment**

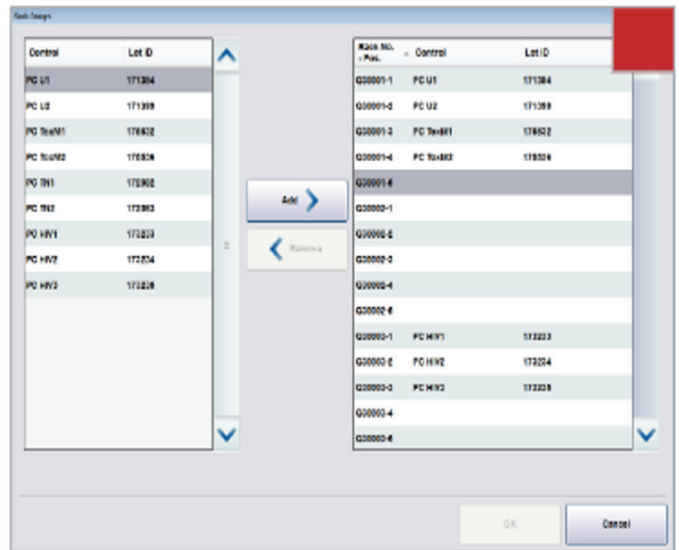


The screenshot shows the 'QC' tab selected in the software interface. Below the navigation tabs, there are buttons for 'Status', 'Run Status', 'Controls', and 'Install'. The 'Controls' button is highlighted. Below these buttons is a table with the following data:

Control	Rack No. - Pos.	S. Type	Lot ID	Exp.
PC Th1		Ser/PI	172082	2015/
PC Th2		Ser/PI	172083	2015/
PC U1	G00001-1	Ser/PI	171394	2015/
PC U2	G00001-2	Ser/PI	171399	2015/
PC Test1	G00001-3	Ser/PI	178532	2015/

Below the table is a large empty area and a 'Rack Assign' button at the bottom right.

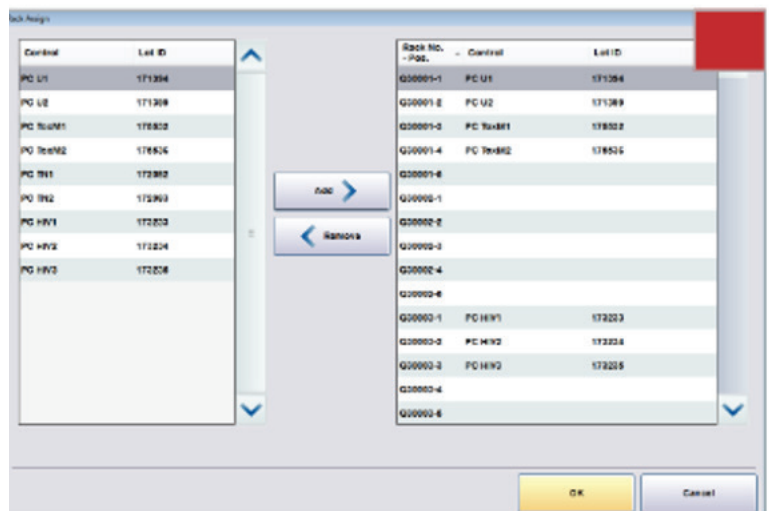
2. On the **Rack Assignment** window select the control from the list on the left. From the right list, choose an unassigned rack ID and rack position. Choose **Add**.



The screenshot shows the 'Rack Assign' dialog box. It has two lists of controls. The left list shows controls with their Lot IDs. The right list shows controls with their Rack No. - Pos. and Lot ID. The 'Add' button is highlighted.

Control	Lot ID
PC U1	171394
PC U2	171399
PC Test1	178532
PC Test2	178536
PC Th1	172082
PC Th2	172083
PC HV1	172203
PC HV2	172204
PC HV3	172205

Rack No. - Pos.	Control	Lot ID
G00001-1	PC U1	171394
G00001-2	PC U2	171399
G00001-3	PC Test1	178532
G00001-4	PC Test2	178536
G00001-6		
G00002-1		
G00002-2		
G00002-4		
G00002-6		
G00003-1	PC HV1	172203
G00003-2	PC HV2	172204
G00003-3	PC HV3	172205
G00003-4		
G00003-6		



The screenshot shows the 'Rack Assign' dialog box with the 'Add' button highlighted. The lists of controls are the same as in the previous screenshot.

Control	Lot ID
PC U1	171394
PC U2	171399
PC Test1	178532
PC Test2	178536
PC Th1	172082
PC Th2	172083
PC HV1	172203
PC HV2	172204
PC HV3	172205

Rack No. - Pos.	Control	Lot ID
G00001-1	PC U1	171394
G00001-2	PC U2	171399
G00001-3	PC Test1	178532
G00001-4	PC Test2	178536
G00001-6		
G00002-1		
G00002-2		
G00002-4		
G00002-6		
G00003-1	PC HV1	172203
G00003-2	PC HV2	172204
G00003-3	PC HV3	172205
G00003-4		
G00003-6		

## Different designation for control racks

There are the following designations for the control racks:



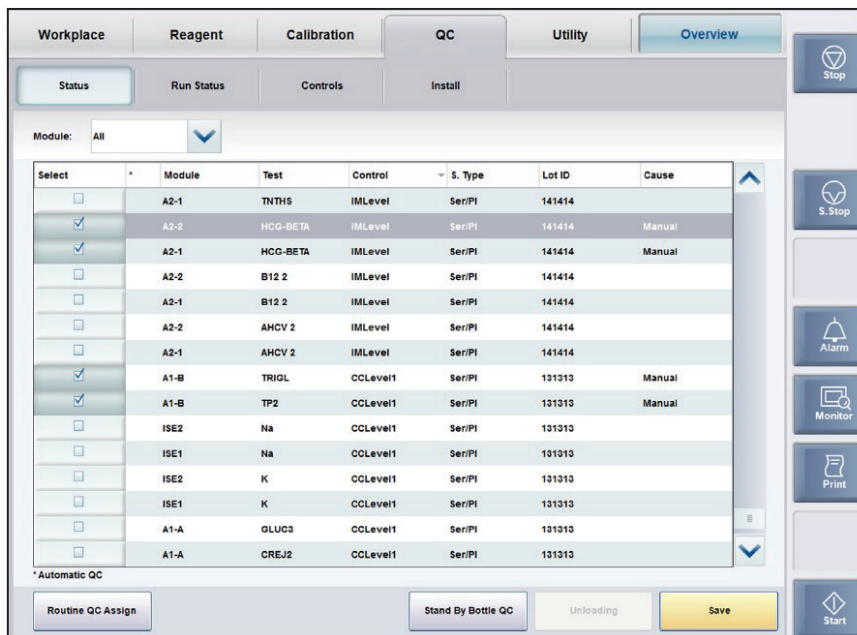
- **Rack Color: White**
- **Label on rack: Q001-Q999**
- **Software displays: Q30001-Q30999**

3. To remove an existing assignment, select the rack number concerned and the rack position from the list on the right and choose **Remove**.
4. Choose **OK** to save the changes and close the window.

## 2. Load

Load IntelliQ QC and run sample

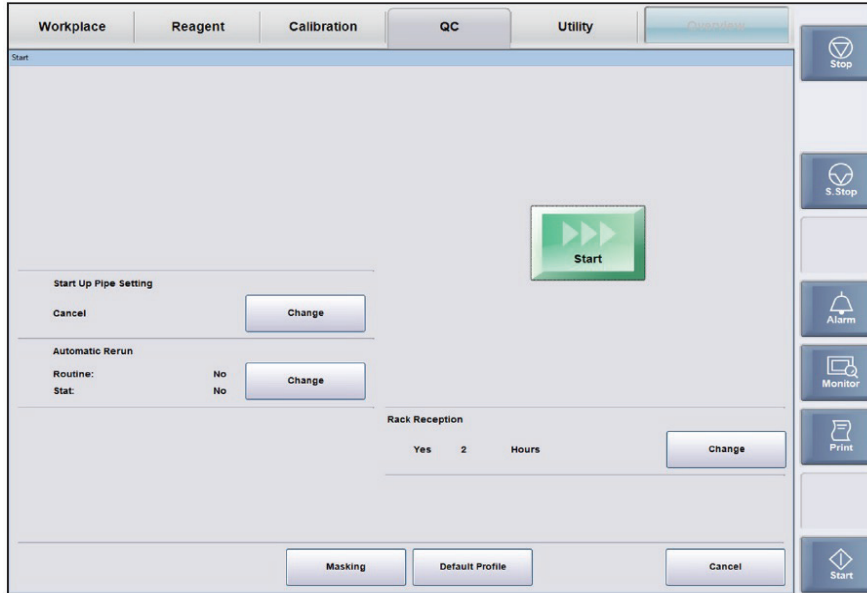
1. Click the **QC tab**.
2. Click the **Status** tab.
3. Highlight desired tests and select them.
4. Click **Save**.



The screenshot shows the software interface with the 'QC' tab selected. The 'Status' sub-tab is active. A table lists various test assignments with columns for Select, Module, Test, Control, S. Type, Lot ID, and Cause. The 'Module' dropdown is set to 'All'. The table contains 16 rows of data, with several rows selected (checked). At the bottom, there are buttons for 'Routine QC Assign', 'Stand By Bottle QC', 'Unloading', 'Save', and 'Start'.

Select	Module	Test	Control	S. Type	Lot ID	Cause
<input type="checkbox"/>	A2-1	TNTH5	IMLevel	Ser/PI	141414	
<input checked="" type="checkbox"/>	A2-2	HCG-BETA	IMLevel	Ser/PI	141414	Manual
<input checked="" type="checkbox"/>	A2-1	HCG-BETA	IMLevel	Ser/PI	141414	Manual
<input type="checkbox"/>	A2-2	B12 2	IMLevel	Ser/PI	141414	
<input type="checkbox"/>	A2-1	B12 2	IMLevel	Ser/PI	141414	
<input type="checkbox"/>	A2-2	AHCV 2	IMLevel	Ser/PI	141414	
<input type="checkbox"/>	A2-1	AHCV 2	IMLevel	Ser/PI	141414	
<input checked="" type="checkbox"/>	A1-B	TRIGL	CCLevel1	Ser/PI	131313	Manual
<input checked="" type="checkbox"/>	A1-B	TP2	CCLevel1	Ser/PI	131313	Manual
<input type="checkbox"/>	ISE2	Na	CCLevel1	Ser/PI	131313	
<input type="checkbox"/>	ISE1	Na	CCLevel1	Ser/PI	131313	
<input type="checkbox"/>	ISE2	K	CCLevel1	Ser/PI	131313	
<input type="checkbox"/>	ISE1	K	CCLevel1	Ser/PI	131313	
<input type="checkbox"/>	A1-A	GLUC3	CCLevel1	Ser/PI	131313	
<input type="checkbox"/>	A1-A	CREJ2	CCLevel1	Ser/PI	131313	

5. Place IntelliQ tube into white QC rack and place on the instrument.
6. Press **Start**.



## Appendix: Auto QC Racks

On Cobas 8000 system, it is possible to keep QC racks in the auto QC compartment in the module sample buffer. The instrument can use auto QC racks for automated QC measurement at any time.

QC samples for auto QC measurements are defined by a range of rack IDs. You specify the range on **Utility>System>Rack Assignment**.

An asterisk(\*) in the second column on the QC> Status submenu indicates which QC materials are used as automatic QC and are loaded onto the instrument.



**Bio-Rad  
Laboratories, Inc.**

*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/diagnostics](http://www.bio-rad.com/diagnostics) **U.S.** 1-800-224-6723 **Australia** +61(2)9914-2800 **Austria** +43-1-877-8901 **Belgium** +32(3)710-53-00 **Brazil** +55(31)3689-6600  
**Canada** 1-514-334-4372 **China** +86 21 6169 8500 **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