

Infinity Lithium Application

Beckman Coulter AU400, AU400e, AU640, AU2700, and AU5400 Parameters

Specific Test Parameters									
General	LIH	ISE	Range						
Test Name:	<input type="text" value="Li"/>	Type:	<input type="text" value="Serum"/>	Operation:	<input type="text" value="Yes"/>				
Sample Volume	<input type="text" value="2"/>	μL	Dilution	<input type="text" value="0"/>	μL	OD Limit			
Pre-Dilution Rate	<input type="text" value="10"/>					Min. OD	<input type="text"/>	Max.O	<input type="text"/>
Reagents Volume	R1 Volume	<input type="text" value="150"/>	μL	Dilution	<input type="text" value="0"/>	μL	Reagent OD limit:		
							First	Low	<input type="text" value="-2.00"/>
							Last	Low	<input type="text" value="-2.00"/>
	R2 Volume	<input type="text" value="0"/>	μL	Dilution	<input type="text" value="0"/>	μL	Dynamic Range Low	<input type="text" value="0.1*"/>	High
							Correlation Factor A	<input type="text" value="1"/>	B
Wavelength:	Pri.	<input type="text" value="520"/>		Sec.	<input type="text" value="480"/>				
Method:	<input type="text" value="END"/>								
Reaction slope:	<input type="text" value="-"/>								
Measuring Point 1:	First	<input type="text" value="0"/>		Last	<input type="text" value="6"/>		Onboard Stability	<input type="text" value="14"/>	Days
Measuring Point 2:	First	<input type="text"/>		Last	<input type="text"/>			<input type="text"/>	
Linearity:	<input type="text"/>								
No Lag Time:	<input type="text"/>								

Specific Test Parameters										
General	LIH	ISE	Range							
Test Name:	<input type="text" value="LI"/>	Type:	<input type="text" value="Serum"/>							
Value/Flag:	<input type="text" value="VALUE"/>	Level L:	<input type="text" value="#"/>			Level H:	<input type="text" value="#"/>			
Specific Ranges:										
	Sex	Year	Month	Year	Month	Low	High			
1.	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>			
2.	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>			
3.	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>			
4.	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>			
5.	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>			
6.	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>			
7.	None selected					<input type="text" value="#"/>	<input type="text" value="#"/>			
8.	Out of Range					<input type="text" value="#"/>	<input type="text" value="#"/>			
Unit	<input type="text" value="mmol/L"/>		Decimal Places	<input type="text" value="#"/>						

Panic Value

Low	High
<input type="text" value="#"/>	<input type="text" value="#"/>

Continued on next page

Infinity Lithium Application

Beckman Coulter AU400, AU400e, AU640, AU2700, and AU5400 Parameters

Calibration Specific					
General		ISE			
Test Name:	<input type="text" value="LI"/>	Type:	<input type="text" value="Serum"/>	<input type="checkbox"/>	
Calibration Type:	<input type="text" value="AB"/>	Formula:	<input type="text" value="Y=AX+B"/>	Counts:	<input type="text" value="#"/>
<Calibrator Parameters>					
	Calibrator	OD	Conc	Factor Range	
	#		†*	Low	High
Point 1:				3.0*	7.0*
Point 2:					
Point 3:					
Point 4:					
Point 5:					
Point 6:					
Point 7:					
1-Point Calibration Point	<input type="text"/>	<input type="checkbox"/>	With CONC-0	Slope Check	<input type="text" value="None"/> ▾
				Advanced Calibration	<input type="text" value="##"/> ▾
MB Type Factor	<input type="text"/>			Calibration Stability	<input type="text" value="7"/> Day

User-defined values
 * To work in SI units (mEq/L) multiply by 1.0

! Do Not Use Lithium Heparin Plasma Samples
 † Beckman Coulter 1.00 mmol/L Calibrator included in kit



WMDE
 Bergerweg 18
 6085 AT Horn
 The Netherlands