

Infinity Lithium Application

Beckman Coulter AU480 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 D	<input type="text"/>
Reagents Volume	R1 (R1-1)	<input type="text" value="150"/>	μL	Dilution	<input type="text" value="0"/>	μL	Reagent OD limit:		
							First Low	<input type="text" value="-2.000"/>	High
							Last Low	<input type="text" value="-2.000"/>	High
	R2 (R2-1)	<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
							Factor for Maker A	<input type="text" value="1"/>	B
Wavelength:	Pri.	<input type="text" value="520"/>		Sec.	<input type="text" value="480"/>		Onboard Stability	<input type="text" value="14"/>	Days
Method:		<input type="text" value="END"/>						<input type="text" value="0"/>	Hour
Reaction slope:		<input type="text" value="-"/>					LIH Influence Check	<input type="text" value="#"/>	
Measuring Point 1:	First	<input type="text" value="0"/>		Last	<input type="text" value="6"/>		Lipemia	<input type="text" value="++++"/>	
Measuring Point 2:	First	<input type="text"/>		Last	<input type="text"/>		Icterus	<input type="text" value="++++"/>	
Linearity:		<input type="text"/>	%				Hemolysis	<input type="text" value="++++"/>	
Lag Time Check:		<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:		From		To		Low	High
	Sex	Year	Month	Year	Month		
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.	No demographics					<input type="text" value="#"/>	<input type="text" value="#"/>
8.	Not within expected values					<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 AU480 Parameters

Calibration Specific										
General		ISE								
Test Name:	<input type="text" value="LI"/>	Type:	<input type="text" value="Serum"/>	<input type="checkbox"/> Use Serum Cal.						
Calibration Type:	<input type="text" value="AB"/>	Formula:	<input type="text" value="Y=AX+B"/>	Counts:	<input type="text" value="#"/>	Slope Check <input type="text" value="None"/>				
<Calibrator Parameters>										
	Calibrator	OD	Conc	Factor Range						
				Low	High					
Point 1:	<input type="text" value="#"/>		<input type="text" value="†*"/>	<input type="text" value="3.0*"/>	<input type="text" value="7.0*"/>					
Point 2:										
Point 3:										
Point 4:										
Point 5:										
Point 6:										
Point 7:										
Point 8:										
Point 9:										
Point 10:										
<Point Cal for Master Curve>										
	Calibrator	OD	Conc	OD Range		No. of Correction Points	<input type="text"/>	Use Master Curve	<input type="text"/>	<input type="checkbox"/> Lot Calibration
				Low	High			Stability		
Point 1:								Reagent Blanks	<input type="text" value="7"/>	Day <input type="text" value="0"/> Hour
Point 2:								Calibration	<input type="text" value="7"/>	Day <input type="text" value="0"/> Hour
MB Type Factor:			1-Point Calibration Point					<input type="checkbox"/> With CONC-0		

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