

# 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 D	<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"/>	High	<input type="text" value="2.500"/>
							Last Low	<input type="text" value="-2.00"/>	High	<input type="text" value="2.500"/>
	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	<input type="text" value="5.0*"/>
Wavelength: Pri.	<input type="text" value="520"/>		Sec.	<input type="text" value="480"/>		Correlation Factor A	<input type="text" value="1"/>	B	<input type="text" value="0"/>	
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 <input type="text"/>					
Measuring Point 2:	First	<input type="text"/>	Last	<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:	<table border="1" style="width: 100%; border-collapse: collapse; text-align: center;"> <thead> <tr> <th></th> <th>Sex</th> <th>Year</th> <th>Month</th> <th>Year</th> <th>Month</th> <th>Low</th> <th>High</th> </tr> </thead> <tbody> <tr><td>1.</td><td>#</td><td>#</td><td>#</td><td>#</td><td>#</td><td>#</td><td>#</td></tr> <tr><td>2.</td><td>#</td><td>#</td><td>#</td><td>#</td><td>#</td><td>#</td><td>#</td></tr> <tr><td>3.</td><td>#</td><td>#</td><td>#</td><td>#</td><td>#</td><td>#</td><td>#</td></tr> <tr><td>4.</td><td>#</td><td>#</td><td>#</td><td>#</td><td>#</td><td>#</td><td>#</td></tr> <tr><td>5.</td><td>#</td><td>#</td><td>#</td><td>#</td><td>#</td><td>#</td><td>#</td></tr> <tr><td>6.</td><td>#</td><td>#</td><td>#</td><td>#</td><td>#</td><td>#</td><td>#</td></tr> <tr><td>7.</td><td colspan="5">None selected</td><td>#</td><td>#</td></tr> <tr><td>8.</td><td colspan="5">Out of Range</td><td>#</td><td>#</td></tr> </tbody> </table>										Sex	Year	Month	Year	Month	Low	High	1.	#	#	#	#	#	#	#	2.	#	#	#	#	#	#	#	3.	#	#	#	#	#	#	#	4.	#	#	#	#	#	#	#	5.	#	#	#	#	#	#	#	6.	#	#	#	#	#	#	#	7.	None selected					#	#	8.	Out of Range					#	#
	Sex	Year	Month	Year	Month	Low	High																																																																										
1.	#	#	#	#	#	#	#																																																																										
2.	#	#	#	#	#	#	#																																																																										
3.	#	#	#	#	#	#	#																																																																										
4.	#	#	#	#	#	#	#																																																																										
5.	#	#	#	#	#	#	#																																																																										
6.	#	#	#	#	#	#	#																																																																										
7.	None selected					#	#																																																																										
8.	Out of Range					#	#																																																																										
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