

Infinity Lithium Application Beckman Coulter AU5800 Parameters

Parameters		Specific Test Parameters					
General	LIH	ISE	Calculated Tests		Range		
Test Name: <input type="text" value="LI"/> ▾		< >		Type: <input type="text" value="Serum"/> ▾	Operation: <input type="text" value="Yes"/> ▾		
Sample Volume	Volume <input type="text" value="2"/> μL	Dilution <input type="text" value="0"/> ▾	OD Limit				
Pre-Dilution Rate <input type="text" value="10"/> ▾				Min OD <input type="text"/>	Max OD <input type="text"/>	<input type="text"/>	
Reagent Volume	R1(R1-1) <input type="text" value="150"/> μL	Dilution <input type="text" value="0"/> μL	Reagent OD Limit				
	R1-2 <input type="text"/>	Dilution <input type="text"/>	First Low <input type="text" value="-2.000"/>	High <input type="text"/>	<input type="text" value="2.500"/>		
	R2(R2-1) <input type="text" value="0"/> μL	Dilution <input type="text" value="0"/> μL	Last Low <input type="text" value="-2.000"/>	High <input type="text"/>	<input type="text" value="2.500"/>		
Common Reagent	Type <input type="text" value="None"/>	Name <input type="text" value="None"/>					
Wavelength:	Pri. <input type="text" value="520"/> ▾nm	Sec. <input type="text" value="480"/> ▾nm					
Method:	<input type="text" value="END"/> ▾						
Reaction slope:	<input type="text" value="-"/> ▾		Dynamic Range	Low <input type="text" value="0.1"/>	High <input type="text"/>	<input type="text" value="5.0"/>	
Measuring Point 1	First <input type="text" value="0"/>	Last <input type="text" value="6"/>	Correlation Factor	A <input type="text" value="1"/>	B <input type="text"/>	<input type="text" value="0"/>	
Measuring Point 2	First <input type="text"/>	Last <input type="text"/>	Factor for Maker	A <input type="text" value="1"/>	B <input type="text"/>	<input type="text" value="0"/>	
Linearity:	<input type="text"/> %						
Lag Time Check	<input type="text"/> ▾		On-board Stability Period	<input type="text" value="14"/> Day	<input type="text" value="#"/> Hour	<input type="text"/>	
LIH Influence Check						<input type="text" value="#"/>	
Lipemia						<input type="text" value="++++"/>	
Icterus						<input type="text" value="++++"/>	
Hemolysis						<input type="text" value="++++"/>	

Parameters		Specific Test Parameters					
General	LIH	ISE	Calculated Tests		Range		
Test Name: <input type="text" value="LI"/> ▾		< >		Type: <input type="text" value="Serum"/> ▾			
Value/Flag: <input type="text" value="#"/> ▾							
Level				Low <input type="text" value="#"/>	High <input type="text" value="#"/>		
Specific Ranges:							
		From		To			
	Sex <input type="text" value="#"/> ▾	Year <input type="text" value="#"/>	Month <input type="text" value="#"/>	Year <input type="text" value="#"/>	Month <input type="text" value="#"/>	Low <input type="text" value="#"/>	
<input type="checkbox"/>	1. <input type="text" value="#"/> ▾	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	High <input type="text" value="#"/>	
<input type="checkbox"/>	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="checkbox"/>	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="checkbox"/>	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="checkbox"/>	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="checkbox"/>	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="#"/>	
7. Standard Demographics				<input type="text"/>	<input type="text"/>		
8. Not within expected values				<input type="text"/>	<input type="text"/>		
Panic Value	Low <input type="text" value="#"/>	High <input type="text" value="#"/>	Unit <input type="text" value="mmol/L"/>		Decimal Places <input type="text" value="#"/>		

Infinity Lithium Application Beckman Coulter AU5800 Parameters

Parameters		Calibration Parameters			
Calibrators	Calibration Specific	STAT Table Calibration			
Test Name: <input type="text" value="LI"/> ▾	<input type="text" value="<"/> <input type="text" value=">"/>	Type: <input type="text" value="Serum"/> ▾	Cuvette <input type="text" value=""/> ▾		
<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="#"/>	
< Calibrator Parameters >					
	Calibrator	OD	Conc	Factor Range Low High	Slope <input type="text" value="None"/> ▾
Point-1	<input type="text" value="#"/> ▾		<input type="text" value="†*"/>	<input type="text" value="3.0"/> <input type="text" value="7.0*"/>	Allowable Range Check <input type="checkbox"/> Reagent Blank <input type="text" value=""/> <input type="checkbox"/> Calibration <input type="text" value=""/> Advanced Calibration Operation <input type="text" value="#"/> ▾ Interval (RB/ACAL) <input type="text" value=""/> ▾ <input type="checkbox"/> Lot Calibration
Point-2					
Point-3					
Point-4					
Point-5					
Point-6					
Point-7					
Point-8					
Point-9					
Point-10					
<Point Cal. For Master Curve>		No. of Correction Points <input type="text" value=""/> ▾	Use Master Curve <input type="text" value=""/> ▾		
	Calibrator	OD	Conc	OD Range Low High	Stability
Point-1	<input type="text" value=""/> ▾				Reagent Blank <input type="text" value="7"/> Day <input type="text" value="#"/> Hour
Point-2	<input type="text" value=""/> ▾				Calibration <input type="text" value="7"/> Day <input type="text" value="#"/> Hour
MB Type Factor <input type="text" value=""/>		1-Point Calibration Point <input type="text" value=""/> ▾		<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