



**Name: Lithium, Dx C 700 AU Lithium Application**

System Reagent: OSRT056  
Test name: LTM1G

Reagent ID: 600

General	LIH	ISE	Calculated Test	Range
<b>Test Name:</b> LTM1G				
<b>Type:</b> Serum				
<b>Operation</b>				
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"/>
Reagent Volume R1 (R1-1)	<input type="text" value="150"/> µL	Dilution	<input type="text" value="0"/> µL	Reagent OD Limit 1 <sup>st</sup> Low <input type="text" value="-2.000"/> High <input type="text" value="2.500"/>
R1-2	<input type="text"/> µL	Dilution	<input type="text"/> µL	Last Low <input type="text" value="-2.000"/> High <input type="text" value="2.500"/>
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 <input type="text" value="5.0*"/>
Common Reagent Type	<input type="text"/>	Name	<input type="text"/>	Correlation Factor A <input type="text" value="1"/> B <input type="text" value="0"/>
Wavelength Pri	<input type="text" value="520"/> nm	Sec	<input type="text" value="480"/> nm	Manufacturer 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="-"/>			Onboard Stability Period <input type="text" value="14"/> Day <input type="text" value="0"/> Hour
Measuring Point-1 1st	<input type="text" value="0"/>	Last	<input type="text" value="6"/>	LIH Influence Check <input type="text" value="#"/>
Measuring Point-2 1st	<input type="text"/>	Last	<input type="text"/>	Lipemia <input type="text" value="£"/>
Linearity Limit	<input type="text"/> %			Icterus <input type="text" value="£"/>
Lag Time Check	<input type="text"/>			Hemolysis <input type="text" value="£"/>

General	LIH	ISE	Calculated Test	Range
<b>Test Name:</b> LTM1G				
<b>Type:</b> Serum				
Value Flag	<input type="text" value="#"/>	Level	Low <input type="text" value="#"/> High <input type="text" value="#"/>	
<b>Specific Ranges</b>				
	Sex	Year	Month	Year
	From	To	Month	Other Type
<input type="checkbox"/> 1:	<input type="text" value="#"/>	<input type="text" value="#"/>	<input type="text" value="#"/>	<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="checkbox"/> 3:	<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="checkbox"/> 5:	<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="#"/>
7:	Standard demographics			<input type="text" value="#"/>
8:	Not within expected values			<input type="text" value="#"/>
Panic Value	Low <input type="text" value="#"/>	High <input type="text" value="#"/>	Unit <input type="text" value="mmol/L*"/>	Select <input type="text" value="Select"/> Decimal Places <input type="text" value="#"/>

Calibrators	General	ISE
<b>Test Name:</b> LTM1G		
<b>Type:</b> 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="#"/>
<Calibrator Parameters>		Slope Check <input type="text" value="None"/>
	Calibrator	OD
		Conc
		Range
		Low
		High
Point-1	<input type="text" value="#"/>	
Point-2	<input type="text"/>	
Point-3	<input type="text"/>	
Point-4	<input type="text"/>	
Point-5	<input type="text"/>	
Point-6	<input type="text"/>	
Point-7	<input type="text"/>	
MB Type Factor	<input type="text"/>	1-Point Calibration Point <input type="text"/>
		<input type="checkbox"/> with Conc-0
		Stability
		Reagent Blank <input type="text" value="7"/> Day <input type="text" value="0"/> Hour
		Calibration <input type="text" value="7"/> Day <input type="text" value="0"/> Hour

# User Defined  
 \* Values set for working in mmol/L. To work in SI units multiply by 1.0  
 C – Closed, Value/Text must be defined in the setting sheet  
 P - Pre-populate, editable, Value/Text must be defined in the setting sheet  
 ±- Sample is only Serum or EDTA plasma.  
 £- Refer to IFU for interference information.  
 ±±- Beckman Coulter 1.00 mmol/L Calibrator included in kit