

Olympus AU2700® / AU5400®

CAT NO:  
TR66056

APPLICATION PROCEDURE

Reagent Preparation: Ready-to-use

Specific Test Parameters			
General	LIH	ISE	Range
Test Name:	LI	<	>
Type:	Serum	Operation: Yes	
Sample Volume:	2 μL	Dil. Volume:	0 μL
Reagent 1 Volume:	150 μL	Dil. Volume:	0 μL
Reagent 2 Volume:	0 μL	Dil. Volume:	0 μL
Wavelength:	Primary: 520	Sec.:	480
Method:	END	Pre-dilution Rate: 10	
Reaction Slope:	NEG	Min OD: L H	
Measuring Point 1:	First: 0	Last: 6	Max OD: L H
Measuring Point 2:	First:	Last:	Reagent OD Limit:
Linearity:	First: %	First: L -2.0000 H 2.5000	
No Lag Time:		Last: L -2.0000 H 2.5000	
		Dynamic Range: L 0.04 H 3.00	
		Correlation Factor: A 1 B 0	
		On Board Stability Period: #	

Specific Test Parameters			
General	LIH	ISE	Range
Test Name:	LI	<	>
Type:	Serum	Value/Flag: #	
		Level L:	#
		Level H:	#
Normal Range:	Age L	Age H	
1 - 6	Sex: #	Year: #	Month: #
7	None selected	Year: #	Month: #
8	Out of range	Year: #	Month: #
Panic Value	L: #	H: #	Unit: mmol/L
		Decimal Places:	2

Calibration Specific			
General	ISE		
Test Name:	LI	<	>
Type:	Serum	Counts: 2	
Cal type:	AB	Formula:	Y = AX + B
Point 1	Cal No. #	OD	Conc
Point 2			
Point 3			
Point 4			
Point 5			
Point 6			
Point 7			
1 - Point cal. Point			With CONC-0
MB type factor		Calibration Stability Period: #	

# User defined parameters

⊛ see recommended LITHIUM Standard TR66901

Olympus AU2700® / AU5400®

**APPLICATION PROCEDURE**

**Reagent Preparation:** Ready-to-use

**CAT NO:**  
TR66056

**\*This is a preliminary instrument application and has not been validated or verified on the Olympus AU2700®/AU5400® analyzer. Each laboratory using this application is responsible for validating and verifying the performance of the application on their system.**

For further information on the use of this reagent, please consult the product insert, or contact Technical Support:  
U.S. 1-800-528-0494, Option 2, International 1-540-869-3200

Fisher Diagnostics, a division of Fisher Scientific, a part of Thermo Fisher Scientific, 8365 Valley Pike, Middletown,  
VA 22645 USA

© 2009 Thermo Fisher Scientific Inc. All rights reserved. AU is a trademark of Olympus America, Inc. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. [www.thermo.com/diagnostics](http://www.thermo.com/diagnostics)