

# DRI™ T-UPTAKE APPLICATION

## Beckman Coulter DxC 700 AU®

Beckman Coulter Reagent REF 0723

The homogeneous enzyme immunoassay is intended to be used for the quantitative determination of unsaturated binding sites on the thyroid binding proteins in human serum or plasma.



For In Vitro Diagnostic Use Only  
Rx Only

### Purpose



The information provided in this application sheet is intended as a supplement to the package insert. Refer to the package insert for information on intended use, reagent storage, reagent preparation, specimen collection, specimen preparation, specimen storage, quality control, and additional performance data. For package inserts, visit [www.thermoscientific.com/diagnostics](http://www.thermoscientific.com/diagnostics) and enter the assay name in the *Search* field.

### Ordering Information

Item	Size	Beckman Coulter Reorder Number
DRI T-Uptake Assay	R1: 1 x 100 mL R2: 1 x 34 mL	0723
DRI T-Uptake Calibrators	5 x 2 mL per level	0738
AU Bottle	20 x 30 mL	63094

### Technical Support

For Technical Support, please contact your local Beckman Coulter Representative.

### Reagent Storage

Refer to the package insert for information on reagent storage. For package inserts, visit [www.thermoscientific.com/diagnostics](http://www.thermoscientific.com/diagnostics) and enter the assay name in the *Search* field.

*Continued on next page*

## Instructions For Use

---

### Procedure for Analyzer

Refer to the operator's manuals for information on analyzer operation. Refer to the package insert for complete reagent preparation.

Prior to pouring into AU bottles, allow the reagent to equilibrate for 15 minutes at refrigerated temperature (2 to 8°C). Dispense R1 reagent and R2 reagent into appropriate AU bottles as shown in the table below.

DRI T-Uptake Assay Kit	AU Reagent Bottle	
	R1 Compartment	R2 Compartment
Enzyme Conjugate Reagent [R1]	One Bottle (30 mL)	
Antibody/Substrate Reagent [R2]		One Bottle (30 mL)

**NOTE:** The Enzyme Conjugate (100 mL kit reagent) is placed in the R1 compartment, and the Antibody/Substrate (34 mL kit reagent) is placed in the R2 compartment. This is the reverse of all other DRI applications.

Warning: These reagents have to be programmed to fixed positions. Do not use the Thermo reagent bottles directly on the AU analyzer.

---

### Results and Data Interpretation

Results for samples will be printed in % T-Uptake.

---

### Specimen Preparation

Refer to the package insert for the complete specimen preparation. The product insert can be found at the Thermo Fisher Scientific website: For package inserts, visit [www.thermoscientific.com/diagnostics](http://www.thermoscientific.com/diagnostics) and enter the assay name in the Search field.

---

### Calibration

Use the DRI T-Uptake Calibrator kit. The calibrators are liquid and ready-to-use.

---

## DRI T-Uptake Assay Beckman Coulter System Parameters, DxC 700 AU

General		LIH		ISE		Calculated Test		Range			
<b>Test Name:</b>		#		Type:		Serum		Operation Yes			
Sample Volume	<input type="text" value="8.0"/>	μL	Dilution	<input type="text" value="0"/>	μL	OD Limit	Min. OD	<input type="text" value="-2.0000"/>	Max OD	<input type="text" value="3.0000"/>	
Pre-Dilution Rate	<input type="text" value="1"/>					Reagent OD Limit	1 <sup>st</sup>	Low	<input type="text" value="-2.0000"/>	High	<input type="text" value="3.0000"/>
Reagent Volume	R1 (R1-1)	<input type="text" value="180"/>	μL	Dilution	<input type="text" value="0"/>	μL	Last	Low	<input type="text" value="-2.0000"/>	High	<input type="text" value="3.0000"/>
	R1-2	<input type="text"/>	μL	Dilution	<input type="text"/>	μL		Low	<input type="text" value="-2.0000"/>	High	<input type="text" value="3.0000"/>
	R2 (R2-1)	<input type="text" value="60"/>	μL	Dilution	<input type="text" value="0"/>	μL	Analytical Measuring Range	Low	<input type="text" value="15.00"/>	High	<input type="text" value="50.00"/>
Common Reagent	Type	<input type="text" value="None"/>		Name	<input type="text" value="None"/>		Correlation Factor	A	<input type="text" value="1"/>	B	<input type="text" value="0"/>
Wavelength	Pri	<input type="text" value="340"/>	nm	Sec	<input type="text" value="520"/>	nm	Manufacturer Factor	A	<input type="text" value="1"/>	B	<input type="text" value="0"/>
Method	<input type="text" value="FIXED1"/>										
Reaction Slope	<input type="text" value="+"/>				Onboard Stability Period		<input type="text" value="30"/>	Day	<input type="text" value="0"/>	Hour	
Measuring Point-1	1st	<input type="text" value="14"/>	Last		<input type="text" value="20"/>	LIH Influence Check		<input type="text" value="No"/>			
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" value="No"/>				Hemolysis		<input type="text" value="+"/>				

General		LIH		ISE		Calculated Test		Range		
<b>Test Name:</b>		#		Type:		Serum				
Value/Flag	<input type="text" value="#"/>		Level		Low	<input type="text" value="#"/>	High	<input type="text" value="#"/>		
Specific Ranges										
	Sex	Year	Month	Year	Month	Other Type	Low	High		
<input type="checkbox"/> 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="None"/>	<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="text" value="#"/>	<input type="text" value="None"/>	<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="None"/>	<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="None"/>	<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="None"/>	<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="None"/>	<input type="text" value="#"/>	<input type="text" value="#"/>		
7:	Standard demographics						<input type="text" value="#"/>	<input type="text" value="#"/>		
8:	Not within expected values						<input type="text" value="#"/>	<input type="text" value="#"/>		
Critical Limits	Low	<input type="text" value="#"/>		High	<input type="text" value="#"/>		Unit	<input type="text" value="%"/>		Select
								Decimal Places	<input type="text" value="#"/>	

# User-defined

## DRI T-Uptake Assay Beckman Coulter System Parameters, DxC 700 AU, *continued*

Calibrators	General	ISE				
<b>Test Name:</b> # <b>Type:</b> Serum						
<input type="checkbox"/> Use Serum Cal.						
Calibration Type: <input type="text" value="5AB"/> Formula: <input type="text" value="EIA Type3"/> Counts: <input type="text" value="2"/>						
<Calibrator Parameters> <span style="float: right;">Slope Check <input type="text" value="+"/></span>						
	Calibrator	OD	Conc	Range		Allowable Range Check <input type="checkbox"/> Reagent Blank <input type="checkbox"/> Calibration Advanced Calibration Operation <input type="text" value="No"/> Interval (RB) <input type="text"/> Interval (ACAL) <input type="text"/>  Stability Reagent Blank <input type="text" value="#"/> Day <input type="text" value="0"/> Hour  Calibration <input type="text" value="#"/> Day <input type="text" value="0"/> Hour
				Low	High	
Point-1	#		15	-2.0000	3.0000	
Point-2	#		20	-2.0000	3.0000	
Point-3	#		30	-2.0000	3.0000	
Point-4	#		40	-2.0000	3.0000	
Point-5	#		50	-2.0000	3.0000	
Point-6						
Point-7						
MB Type Factor <input type="text"/> 1-Point Calibration Point <input type="text" value="None"/> <input type="checkbox"/> with Conc-0						

# User-defined

## Additional Information

### Important

Since Beckman Coulter does not manufacture the reagent or perform quality control or other tests on individual lots, Beckman Coulter cannot be responsible for the quality of the data obtained which is caused by performance of the reagent, any variation between lots of reagent, or protocol changes by the Manufacturer.

### Shipping Damage

Please notify your Beckman Coulter Technical Support Center if this product is received damaged.

© 2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. AU Series Systems are trademarks of Beckman Coulter. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others.

*End*