

DRI™ T-UP TAKE APPLICATION

Beckman Coulter DxC 500 AU / 500i

Beckman Coulter Reagent REF 0723

The Application is Intended 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

Intended Use



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.thermofisher.com 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 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 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. Refer to the package insert for the concentration of each calibrator.

Reagent Name: DRI T-Uptake Assay REF 0723 DxC 500 AU / 500i Serum (Plasma)
Settings
Calibrator Name: DRI T-Uptake Calibrator Kit REF 0738

Reagent ID 564

TEST CONFIGURATION & CHEMISTRY DETAILS

Assay Name	Test	Rev	Discipline	Chemistry		
Test ID	cTUP		Calculated Result	<input type="checkbox"/>		
LIS Code	cTUP		Result Type	Quantitative ▼		
UNITS AND RANGE SETTINGS						
Use Settings from	Serum ▼	Units	% ▼	Decimal Places	x.xx ▼	Plasma
Test Kind	General ▼	Revision	01	<input checked="" type="checkbox"/> Multi Reagent Switch		
Reagent Name	cTUP	Reagent ID	564	<input type="checkbox"/> FSE Test		
ABB Name	T-U1G	Parameter Long Name	T-Uptake 564 T-U1G Serum			
Region	<input checked="" type="checkbox"/> US	<input checked="" type="checkbox"/> OUS	<input checked="" type="checkbox"/> AP	<input type="checkbox"/> JP	<input checked="" type="checkbox"/> EU	<input type="checkbox"/> Other

GENERAL PARAMETERS

SAMPLE VOLUME	Sample Volume	8.0	μL	Dilution	0	μL	REACTION OD LIMIT	Low	-2.0000	High	3.0000		
	Predilution Rate	1					REACTION BLANK OD LIMIT	First: Low	-2.0000	High	3.0000		
REAGENT VOLUME	R1-1	180	μL	Dilution	0	μL		Last: Low	-2.0000	High	3.0000		
	R2-1	60	μL	Dilution	0	μL	ANALYTICAL MEASURING RANGE	Low	15.00	High	50.00		
WAVELENGTH	Primary	340	nm	Secondary	520	nm	MANUFACTURER FACTOR	A	1	B	0		
METHOD	FIXED 1 ▼						REAGENT ONBOARD STABILITY	29		Days	0	Hours	
REACTION SLOPE	+						LIH INFLUENCE CHECK	<input type="checkbox"/> Perform LIH check					
MEASURING POINT	Point 1: First	14	Last	20									
	Point 2: First		Last		Lipemia	+	▼	Icterus	+	▼	Hemolysis	+	▼
Linearity Limit			%										
Lag Time Check			<input type="checkbox"/> Perform Lag Time Check										

Reagent Name: DRI T-Uptake Assay REF 0723 DxC 500 AU / 500i Serum (Plasma)

Reagent ID 564

Settings

Calibrator Name: DRI T-Uptake Calibrator Kit REF 0738, *Continued*

CALIBRATION PARAMETERS

Base Unit	Decimal Place	Unit 1	Factor 1	Unit 2	Factor 2	Unit 3	Factor 3	Unit 4	Factor 4
%	▼ 2	▼ None	▼ 0	None	▼ 0	None	▼ 0	None	▼ 0

CALIBRATOR SPECIFIC

Calibration Type

Counts

CALIBRATION OD AND CONCENTRATION PARAMETERS

☐ Use highest calibrator for Upper AMR

Formula

MB Factor

Calibrator Name

Positive Cutoff

Add

☒ SLOPE CHECK

Number of Levels

Slope Check

STABILITY AND INTERVAL

Reagent Blank Stability Days Hours

Interval

Calibration Stability Days Hours

Interval

OD DELTA CHECK

☐ Reagent Blank

☐ Calibration

	Calibrator Name	Conc	OD Range Low	OD Range High
Point 1	cTUP CAL-1	15.00	-2.00	3.00
Point 2	cTUP CAL-2	20.00	-2.00	3.00
Point 3	cTUP CAL-3	30.00	-2.00	3.00
Point 4	cTUP CAL-4	40.00	-2.00	3.00
Point 5	cTUP CAL-5	50.00	-2.00	3.00
Point 6				
Point 7				

PROZONE CHECK PARAMETERS

☐ Logic Check 1

Check Points

Point 1	<input type="text" value="0"/>
Point 2	<input type="text" value="0"/>
Point 3	<input type="text" value="0"/>

Decision Values

Value 1	<input type="text" value="0"/>
Value 2	<input type="text" value="0"/>
Value 3	<input type="text" value="0"/>

☐ Logic Check 2

Check Points

Point 1	<input type="text" value="0"/>
Interval	<input type="text" value="1"/>

Decision Values

Value 1	<input type="text" value="0"/>
Value 2	<input type="text" value="0"/>

☐ Logic Check 3

Check Points

Point 1	<input type="text" value="0"/>
Interval	<input type="text" value="1"/>

Decision Values

Value 1	<input type="text" value="0"/>
Value 2	<input type="text" value="0"/>

Limit Points

Limit 1	<input type="text" value="0"/>
Limit 2	<input type="text" value="27"/>

Limit Points

Limit 1	<input type="text" value="0"/>
Limit 2	<input type="text" value="27"/>

Limit Points

Limit 1	<input type="text" value="0"/>
Limit 2	<input type="text" value="27"/>

Check Pattern

Pattern

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.

© 2024 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