

QMS™ EVEROLIMUS APPLICATION
Beckman Coulter DxC 700 AU®



Beckman Coulter Reagent REF A53729 (US), A53716 (OUS)

The QMS Everolimus Assay is to be used for the quantitative determination of everolimus in human whole blood on automated clinical chemistry analyzers.



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 and enter the assay name in the *Search* field.

Ordering Information

| Item | Size | Beckman Coulter Reorder Number |
|-------------------------------|--|--------------------------------|
| QMS Everolimus Assay | R1: 1 x 22 mL R2: 1 x 8 mL Precipitation Reagent: 1 x 8 mL | A53729 (US) A53716 (OUS) |
| QMS Everolimus Calibrator Set | 6 levels 1 x 3 mL per level | A53721 (US) A53724 (OUS) |
| QMS Everolimus Control Set | 3 levels 1 x 3 mL per level | A53717 |
| AU Bottle | 20 x 15 mL | 63165 |
| 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:

| QMS Everolimus Assay Kit | AU Reagent Bottle | |
|---|--------------------|--------------------|
| | R1 Compartment | R2 Compartment |
| Anti-Everolimus Polyclonal Antibody R1 | One Bottle (30 mL) | |
| Everolimus-Coated Microparticles R2 | | One Bottle (15 mL) |

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

Significant interference from QMS Everolimus into the Urine/CSF Albumin (B38858/B46435) assays has been observed due to reagent carryover in random access analyzers. Setup the recommended contamination parameters below:

| DxC 700 AU Contamination Parameters | | | | | | | |
|-------------------------------------|------------------------------------|-------------------------|---------------|------------|--------------------|-------------|---------|
| No. | Providing Test Name | Receiving Test Name | Reagent Probe | Wash Count | Effective of Water | Prevent Use | |
| | | | | | | Mixer | Cuvette |
| 1 | EVR (A53729 US) (A53716 OUS) | UALB (B38858/B46435) | Water | 3 | Yes | Yes | Yes |

Results and Data Interpretation

Results for samples will be printed in ng/mL.

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 QMS Everolimus Calibrator set. The calibrators are prepared like patient samples.

QMS Everolimus Assay Beckman Coulter System Parameters, DxC 700 AU

| General | LIH | ISE | Calculated Test | Range |
|---|--|---|--|-------|
| Test Name: # <input type="text"/> Type: Serum Operation: Yes | | | | |
| Sample Volume | <input type="text" value="10"/> μL | Dilution <input type="text" value="0"/> μL | OD Limit | |
| Pre-Dilution Rate | <input type="text" value="1"/> | | Min. OD <input type="text" value="-2.0000"/> Max OD <input type="text" value="3.0000"/> | |
| Reagent Volume | R1 (R1-1) <input type="text" value="175"/> μL | Dilution <input type="text" value="0"/> μL | Reagent OD Limit 1 st 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 | Last Low <input type="text" value="-2.0000"/> High <input type="text" value="3.0000"/> | |
| | R2 (R2-1) <input type="text" value="45"/> μL | Dilution <input type="text" value="0"/> μL | Analytical Measuring Range Low <input type="text" value="1.500"/> High <input type="text" value="20.00"/> | |
| 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="700"/> nm | Sec <input type="text" value="None"/> 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="29"/> Day <input type="text" value="0"/> Hour | |
| Measuring Point-1 | 1st <input type="text" value="24"/> Last <input type="text" value="27"/> | | 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 |
|---|---|--|--|---|
| Test Name: # <input type="text"/> Type: Serum | | | | |
| Value/Flag | <input type="text" value="#"/> | Level | Low <input type="text" value="#"/> High <input type="text" value="#"/> | |
| Specific Ranges | | | | |
| | From | To | Other Type | Low High |
| <input type="checkbox"/> 1: | 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="#"/> | <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="ng/mL"/> | Select <input type="text" value="2"/> Decimal Places |

User-defined

QMS Everolimus Assay Beckman Coulter System Parameters, DxC 700 AU, *continued*

| Calibrators | General | ISE | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | |
|---|------------|-----|-----|------------|--------|------|-------|-------|-----|------|---------|-----|--|---|---------|--------|---------|-----|--|-----|---------|--------|---------|-----|--|---|---------|--------|---------|-----|--|---|---------|--------|---------|-----|--|----|---------|--------|---------|-----|--|----|---------|--------|---------|-----|--|--|--|--|
| Test Name: # ▼ Type: Serum ▼ <input type="checkbox"/> Use Serum Cal. | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | |
| Calibration Type: 6AB ▼ Formula: EIA Type 1 ▼ Counts: 2 | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | |
| <Calibrator Parameters> <table border="1" style="width: 100%;"> <thead> <tr> <th rowspan="2"></th> <th rowspan="2">Calibrator</th> <th rowspan="2">OD</th> <th rowspan="2">Conc</th> <th colspan="2">Range</th> </tr> <tr> <th>Low</th> <th>High</th> </tr> </thead> <tbody> <tr> <td>Point-1</td> <td># ▼</td> <td></td> <td>0</td> <td>-2.0000</td> <td>3.0000</td> </tr> <tr> <td>Point-2</td> <td># ▼</td> <td></td> <td>1.5</td> <td>-2.0000</td> <td>3.0000</td> </tr> <tr> <td>Point-3</td> <td># ▼</td> <td></td> <td>3</td> <td>-2.0000</td> <td>3.0000</td> </tr> <tr> <td>Point-4</td> <td># ▼</td> <td></td> <td>6</td> <td>-2.0000</td> <td>3.0000</td> </tr> <tr> <td>Point-5</td> <td># ▼</td> <td></td> <td>12</td> <td>-2.0000</td> <td>3.0000</td> </tr> <tr> <td>Point-6</td> <td># ▼</td> <td></td> <td>20</td> <td>-2.0000</td> <td>3.0000</td> </tr> <tr> <td>Point-7</td> <td># ▼</td> <td></td> <td></td> <td></td> <td></td> </tr> </tbody> </table> | | | | Calibrator | OD | Conc | Range | | Low | High | Point-1 | # ▼ | | 0 | -2.0000 | 3.0000 | Point-2 | # ▼ | | 1.5 | -2.0000 | 3.0000 | Point-3 | # ▼ | | 3 | -2.0000 | 3.0000 | Point-4 | # ▼ | | 6 | -2.0000 | 3.0000 | Point-5 | # ▼ | | 12 | -2.0000 | 3.0000 | Point-6 | # ▼ | | 20 | -2.0000 | 3.0000 | Point-7 | # ▼ | | | | |
| | Calibrator | OD | | | | | Conc | Range | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | |
| | | | Low | High | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | |
| Point-1 | # ▼ | | 0 | -2.0000 | 3.0000 | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | |
| Point-2 | # ▼ | | 1.5 | -2.0000 | 3.0000 | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | |
| Point-3 | # ▼ | | 3 | -2.0000 | 3.0000 | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | |
| Point-4 | # ▼ | | 6 | -2.0000 | 3.0000 | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | |
| Point-5 | # ▼ | | 12 | -2.0000 | 3.0000 | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | |
| Point-6 | # ▼ | | 20 | -2.0000 | 3.0000 | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | |
| Point-7 | # ▼ | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | |
| MB Type Factor: <input type="text"/> 1-Point Calibration Point: None ▼ <input type="checkbox"/> with Conc=0 | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | |
| Slope Check: - ▼ Allowable Range Check <input type="checkbox"/> Reagent Blank <input type="text"/> <input type="checkbox"/> Calibration <input type="text"/> Advanced Calibration Operation: No ▼ Interval (RB): <input type="text"/> ▼ Interval (ACAL): <input type="text"/> ▼ | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | |
| Stability Reagent Blank: # Day 0 Hour Calibration: # Day 0 Hour | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | | |

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.



EC REP B-R-A-H-M-S GmbH, Neuendorfstrasse 25, 16761, Hennigsdorf, Germany

End