

Olympus AU2700® / AU5400®

CAT NO:  
TR25421

APPLICATION PROCEDURE

Reagent Preparation: Ready-to-use

| Specific Test Parameters |               |                   |                              |
|--------------------------|---------------|-------------------|------------------------------|
| General                  | LIH           | ISE               | Range                        |
| Test Name:               | AMY ▼         | < >               | Type: Serum                  |
|                          |               |                   | Operation: Yes ▼             |
| Sample Volume:           | 3 μL          | Dil. Volume: 0 μL | Pre-dilution Rate: 1         |
| Reagent 1 Volume:        | 120 μL        | Dil. Volume: 0 μL | Min OD: L 0.000 H 2.5000     |
| Reagent 2 Volume:        | μL            | Dil. Volume: 0 μL | Max OD: H 2.5000             |
| Wavelength:              | Primary 410 ▼ | Sec. 520 ▼        | Reagent OD Limit:            |
| Method:                  | RATE          |                   | First: L 0.0000 H 0.7000     |
| Reaction Slope:          | + ▼           |                   | Last: L 0.0000 H 0.7000      |
| Measuring Point 1:       | First 4       | Last 10           | Dynamic Range:               |
| Measuring Point 2:       | First         | Last              | L 10 H 2000                  |
| Linearity:               | First %       |                   | Correlation Factor:          |
| No Lag Time:             | YES ▼         |                   | A 1 B 0                      |
|                          |               |                   | On Board Stability Period: # |

| Specific Test Parameters |               |            |                   |
|--------------------------|---------------|------------|-------------------|
| General                  | LIH           | ISE        | Range             |
| Test Name:               | AMY ▼         | < >        | 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: U/L         |
|                          |               |            | Decimal Places: 0 |

| Calibration Specific |         |                       |                               |
|----------------------|---------|-----------------------|-------------------------------|
| General              | ISE     |                       |                               |
| Test Name:           | AMY ▼   | < >                   | Type: Serum                   |
| Cal type             | MB ▼    | Formula: Y = AX + B ▼ | Counts: 2                     |
|                      | Cal No. | OD                    | Process: # ▼                  |
| Point 1              |         | Conc                  | Factor/OD-L                   |
| Point 2              |         |                       | Factor/OD-H                   |
| Point 3              |         |                       |                               |
| Point 4              |         |                       |                               |
| Point 5              |         |                       |                               |
| Point 6              |         |                       |                               |
| Point 7              |         |                       |                               |
| 1 - Point cal. Point |         | With CONC-0           |                               |
| MB type factor       | 5313    |                       | Calibration Stability Period: |

# User defined parameters

Olympus AU2700® / AU5400®

**APPLICATION PROCEDURE**

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

**CAT NO:**  
TR25421

**\*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)