



CEDIA[®] CYCLOSPORINE PLUS HIGH RANGE APPLICATION

ORTHO CLINICAL DIAGNOSTICS VITROS[®] XT 7600 INTEGRATED SYSTEM, VITROS[®] 5600 INTEGRATED SYSTEM, VITROS[®] 4600 CHEMISTRY SYSTEM, AND VITROS[®] 5,1 FS CHEMISTRY SYSTEM

Reference No. 100147

This Application is Intended for the Quantitative Determination of Cyclosporine in Human Whole Blood



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.orthoclinicaldiagnostics.com > Ortho Care > Technical Documents > MicroTip Partnership Assays (MPA).

Ordering Information

Please place your order with Ortho Clinical Diagnostics. Ordering information available on www.orthoclinicaldiagnostics.com.

Technical Support Information

Contact Ortho Clinical Diagnostics for technical support. Contact information available on www.orthoclinicaldiagnostics.com.



Microgenics Corporation, part of Thermo Fisher Scientific
46500 Kato Road, Fremont, CA 94538 USA
U.S. Toll free: (800) 232-3342 / Tel: (510) 979-5000
U.S. Toll free fax: (888) 527-8001 / Fax: (510) 979-5420



B·R·A·H·M·S GmbH, Neuendorfstrasse 25, 16761 Hennigsdorf, Germany

Reagent Pack Storage

Reconstituted reagents are stable for 60 days at 2-8°C when stored in their original container or in UDxx reagent packs.

Reconstituted reagents stored in UDxx reagent packs onboard the analyzer are stable for 28 days. Reagent life can be extended by storing the reagent packs in a refrigerator at 2-8°C between use.

To split the reconstituted reagent among reagent packs of the same UDxx ID, follow the chart below:

| Number of packs | EA (mL) in UDxx/A | ED (mL) in UDxx/B | Tests/pack |
|-----------------|-------------------|-------------------|------------|
| 3 | 13.7 | 6.3 | 74 |

NOTE: Once the individual UDxx pack number is selected for use during the protocol programming, it is the only UDxx pack number to use for this protocol. The same pack can be used for both the Low Range and High Range Cyclosporine PLUS assays at the same time (each is calibrated separately).

Special Reagent Packs for User Defined Assays

(Please order from Ortho Clinical Diagnostics; not available from Microgenics)

| Part Number | Description | Quantity |
|-------------|---------------------|------------|
| 680 2246 | UD01 Packs (Empty) | 1 BOX/6PKS |
| 680 2247 | UD02 Packs (Empty) | 1 BOX/6PKS |
| 680 2248 | UD03 Packs (Empty) | 1 BOX/6PKS |
| 680 2249 | UD04 Packs (Empty) | 1 BOX/6PKS |
| 680 2250 | UD05 Packs (Empty) | 1 BOX/6PKS |
| 680 2251 | UD06 Packs (Empty) | 1 BOX/6PKS |
| 680 2252 | UD07 Packs (Empty) | 1 BOX/6PKS |
| 680 2253 | UD08 Packs (Empty) | 1 BOX/6PKS |
| 680 2254 | UD09 Packs (Empty) | 1 BOX/6PKS |
| 680 2255 | UD10 Packs (Empty) | 1 BOX/6PKS |
| 684 4449 | UD11 Packs (Empty) | 1 BOX/6PKS |
| 684 4448 | UD12 Packs (Empty) | 1 BOX/6PKS |
| 684 4445 | UD13 Packs (Empty) | 1 BOX/6PKS |
| 684 4442 | UD14 Packs (Empty) | 1 BOX/6PKS |
| 684 4447 | UD15 Packs (Empty) | 1 BOX/6PKS |
| 684 4444 | UD16 Packs (Empty) | 1 BOX/6PKS |
| 684 4441 | UD17 Packs (Empty) | 1 BOX/6PKS |
| 684 4446 | UD18 Packs (Empty) | 1 BOX/6PKS |
| 684 4443 | UD19 Packs (Empty) | 1 BOX/6PKS |
| 684 4440 | UD20 Packs (Empty) | 1 BOX/6PKS |
| 680 2256 | UDDL1 Packs (Empty) | 1 BOX/6PKS |
| 680 2257 | UDDL2 Packs (Empty) | 1 BOX/6PKS |

Calibration Frequency

It is recommended that recalibration occur after reagent pack change, after calibrator lot change, after performance of monthly instrument maintenance and as required following quality control procedure.

Sample Preparation

Follow the CEDIA Cyclosporine PLUS Assay Sample Preparation as described in the package insert.

CEDIA Cyclosporine PLUS High Range Assay
Ortho Clinical Diagnostics VITROS® XT 7600 System, VITROS® 5600 System,
VITROS® 4600 System, and VITROS® 5,1 FS System Parameters

Configure Assay

Full Assay Name: Cyclosporine HR Version Date: 9/15/2010
 Short Assay Name: CsAHR Fluid Type: Wh Blood
 Assay Model Type: 2 Point Rate Template: *2Pt R1-S-R2
 Cal Model Type: Linear Calibrator Bottles: 2 Reagent Reps per Cal : 2

Reagent Lot Information

On-Board Stability: 28 Days
 Reagent Lot Num. Kit Lot
 Shelf Exp. Date: Kit Exp Date

Edit Dilution Parameters

Diluent: None Standard Dilution Factor: 1.0
 Reflex Dilution: OFF Dilution Factor: 1.0
 Reduction Factor: 1.0

Edit Result Parameters

Units: ng/mL Reference Interval: 0.0 to 90000000
 Significant Digits: 4 Precision Digits: 3 Supplementary: 0.0 to 90000000
 User Adjusted Parameters Reportable Range: 450 to 2000
 Slope: 1.0 Intercept: 0.0 **(More Assay Parm) – Edit 2 Pt Rate Additional Parameters**
 CuveTip Exp Time: 35 Temp Sens : No Initial Abs. Limits: -0.20 to 3.50 (2.7 for 5,1 FS*)
 Second Abs. Limits: -0.20 to 3.50 (2.7 for 5,1 FS*)
 Antigen Excess Factor: 9.0

Edit Protocol Parameters

| Step | Volume | Pack ID | Seconds | Wavelength |
|---------------|---------|---------|---------|------------|
| 1. Reagent | 150 uL | UDxx /A | | |
| 2. Incubation | | | 0.0 | |
| 3. Sample | 2.6 uL | | | |
| 4. Incubation | | | 304.0 | |
| 5. Reagent | 66.4 uL | UDxx /B | | |
| 6. Incubation | | | 228.0 | |
| 7. Read | | | | 575 nm |
| 8. Incubation | | | 76.0 | |
| 9. Read | | | | 575 nm |

**CEDIA Cyclosporine PLUS High Range Assay
 Ortho Clinical Diagnostics VITROS® XT 7600 System, VITROS® 5600 System,
 VITROS® 4600 System, and VITROS® 5,1 FS System Parameters, *continued***

Edit Calibration Parameters

| Bottle # | Dil Factor | Cal Rep Resp Range | Calibrator Lot: <u>Cal Kit lot</u> |
|----------|------------|--------------------|------------------------------------|
| 1 | <u>1.0</u> | <u>0.20</u> | Cal value: <u>per cal lot</u> |
| 2 | <u>1.0</u> | <u>0.20</u> | Cal Value: <u>per cal lot</u> |

(More Cal Parm) – Edit Linear or Logit/Log Additional Parameters

| | | |
|-------------------------------|------------------------------|--------------------------------------|
| Monotonicity: <u>Increase</u> | | |
| Max Resp High: <u>3.00</u> | Min. Resp. High: <u>3.00</u> | Cal Fit Goodness Limit: <u>0.990</u> |
| Max Resp. Low: <u>-3.00</u> | Min Resp. Low: <u>-3.00</u> | |

Edit Triple Read Parameters

| | Reportable Conc. | Triple Read Limit |
|------------------|------------------|-------------------|
| Reportable Min.: | <u>450</u> | <u>10</u> |
| Critical Conc.: | <u>1000</u> | <u>8.0</u> % |
| Reportable Max.: | <u>2000</u> | <u>8.0</u> % |

Comments: Please see important assay and analyzer notes on page 2.

*** For the 5,1 FS the upper limit for initial and secondary abs limits is 2.70. Samples with high hematocrit levels may return a cuvette blank error. In these cases, please dilute the sample as per the package insert instructions and re-run.**

© 2019 Thermo Fisher Scientific, Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. VITROS is a registered trademark of Ortho Clinical Diagnostics.

End