

**CEDIA[®] CYCLOSPORINE PLUS LOW RANGE
APPLICATION
ORTHO CLINICAL DIAGNOSTICS VITROS[®] XT 7600
INTEGRATED SYSTEM, VITROS[®] 5600 INTEGRATED
SYSTEM, AND VITROS[®] 4600 AND 5,1 FS CHEMISTRY
SYSTEMS**

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 Low 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 LR Version Date: 24 September 2015
 Short Assay Name: CsALR 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: 25 to 450
 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	16.8 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 Low 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>25</u>	<u>10</u>
Critical Conc.:	<u>200</u>	<u>8.0</u> %
Reportable Max.:	<u>450</u>	<u>8.0</u> %

*** 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