

**DRI™ PROPOXYPHENE APPLICATION**



**ORTHO CLINICAL DIAGNOSTICS VITROS® XT 7600  
INTEGRATED SYSTEM, VITROS® 5600 INTEGRATED  
SYSTEM, AND VITROS® 4600 CHEMISTRY SYSTEM**

Reference No. 0432

Intended for Qualitative and Semiquantitative Determination of Propoxyphene in Human Urine



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](http://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](http://www.orthoclinicaldiagnostics.com).

**Technical  
Support  
Information**

Contact Ortho Clinical Diagnostics for technical support. Contact information available on [www.orthoclinicaldiagnostics.com](http://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

Unopened reagents are stable until the expiration date when stored at 2 to 8°C.

Reagents stored in UDxx reagent packs onboard the analyzer are stable for 30 days.

Reagents are supplied liquid ready-to-use.

The UDxx reagent pack may be filled up to 15 mL of Reagent A in UDxx/A bottle and 15 mL of Reagent B in UDxx/B bottle, and provides 118 tests.

R1 (mL) in UDxx/A	R2 (mL) in UDxx/B	Tests/pack
15	15	118

**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.

### 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

## Results and Data Interpretation

For the Qualitative assay, the VITROS<sup>®</sup> XT 7600 System, VITROS<sup>®</sup> 5600 System, and VITROS<sup>®</sup> 4600 System can report both a unit-less Index number and a NEG / POS classification. The assay is calibrated at the cutoff concentration so that the Index number less than the cutoff value are negative and those greater than the cutoff value are positive.

**NOTE:** Qualitative results must only be reported as Negative or Positive. Unit-less Index numbers cannot be used in reportable results.

For the Semiquantitative assay, the VITROS<sup>®</sup> XT 7600 System, VITROS<sup>®</sup> 5600 System, and VITROS<sup>®</sup> 4600 System can report both Semiquantitative numbers in ng/mL and a NEG/POS classification. The POS classification will also be given on samples above the Semiquantitative reporting range. Sample results above the high calibrator should be diluted with negative urine and retested.

---

## Out of Range Error Codes

A very high analyte sample can produce an absorbance outside the VITROS<sup>®</sup> System photometer range and will result in a CB error code. Sample results above the high calibrator should be diluted with negative urine and retested.

With the Semiquantitative assay, a sample above the logit-log4 spline prediction range, but within the System photometer range, will produce an ER error. Sample results above the high calibrator should be diluted with negative urine and retested.

---

## Calibration Frequency

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

# DRI Propoxyphene Assay – Qualitative Ortho Clinical Diagnostics VITROS® XT 7600 System, VITROS® 5600 System, and VITROS® 4600 System Parameters

Full Assay Name: Propoxy Qual Version Date: 09/05/2017  
Short Assay Name: PP300 Fluid Type: Urine  
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: 30 Days

Reagent Lot Num: Kit Lot

Shelf Exp. Date: Kit Exp Date

## Edit Dilution Parameters

Diluent: DATDIL Standard Dilution Factor: 1.0

Reflex Dilution: OFF Dilution Factor: 1.0

Reduction Factor: 1.0

---

## Edit Result Parameters

**NOTE: For XT 7600 / 5600 / 4600 only enter qualitative ranges and the Measuring (Reportable) range**

Units: \_\_\_\_\_

Number of Qualitative Ranges: 2

Significant Digits: 5 Precision Digits: 1

1. NEGATIVE

User Adjusted Parameters

2. POSITIVE 300 or greater

Slope: 1.0 Intercept: 0.0

Report Qualitative Result Outside of Range: Yes

CuveTip Exp Time: 35 Temp Sens : No

Measuring (Reportable) 0 to 10000

## (More Assay Parm) – Edit 2 Pt Rate Additional Parameters

Initial Abs. Limits: -0.20 to 3.50

Second Abs. Limits: -0.20 to 3.50

Antigen Excess Factor: 9.0

---

**DRI Propoxyphene Assay – Qualitative**  
**Ortho Clinical Diagnostics VITROS® XT 7600 System, VITROS® 5600 System, and**  
**VITROS® 4600 System Parameters, *continued***

**Edit Protocol Parameters**

Step	Volume	Pack ID	Seconds	Wavelength
1. Reagent	100 µL	UDxx /A		
2. Incubation			0.0	
3. Sample	8			
4. Incubation			304.0	
5. Reagent	100 µL	UDxx /B		
6. Incubation			38.0	
7. Read				340 nm
8. Incubation			76.0	
9. Read				340 nm

**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>0</u>
2	<u>1.0</u>	<u>0.20</u>	Cal Value: <u>300</u>

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

Monotonicity: Increase

Max Resp High: 3.00      Min. Resp. High: 3.00      Cal Fit Goodness Limit: 0.990

Max Resp. Low: -3.00      Min Resp. Low: -3.00

**Edit Triple Read Parameters**

	Reportable Conc.	Triple Read Limit
Reportable Min.:	<u>0</u>	<u>24.00</u>
Critical Conc.:	<u>300</u>	<u>8.0</u> %
Reportable Max.:	<u>10000</u>	<u>8.0</u> %

**Comments: N/A**

**DRI Propoxyphene Assay – Semiquantitative**  
**Ortho Clinical Diagnostics VITROS® XT 7600 System, VITROS® 5600 System, and**  
**VITROS® 4600 System Parameters**

Full Assay Name: Propoxy SemiQuant Version Date: 09/05/2017  
Short Assay Name: PPX Fluid Type: Urine  
Assay Model Type: 2 Point Rate Template: \*2Pt R1-S-R2  
Cal Model Type: Logit/Log4 Calibrator Bottles: 5 Reagent Reps per Cal: 2

---

**Reagent Lot Information**

On-Board Stability: 30 Days  
Reagent Lot Num: Kit Lot  
Shelf Exp. Date: Kit Exp Date

**Edit Dilution Parameters**

Diluent: DATDIL Standard Dilution Factor: 1.0  
Reflex Dilution: OFF Dilution Factor: 1.0  
Reduction Factor: 1.0

---

**Edit Result Parameters**

**NOTE: For XT 7600 / 5600 / 4600 only enter qualitative ranges and the Measuring (Reportable) range**

Units: ng/mL Number of Qualitative Ranges: 2  
Significant Digits: 5 Precision Digits: 1  
User Adjusted Parameters  
Slope: 1.0 Intercept: 0.0  
Cuvette Exp Time: 35 Temp Sens : No

1. NEGATIVE  
2. POSITIVE 300 or greater  
Report Qualitative Result Outside of Range: Yes  
Measuring (Reportable) 0 to 1000

**(More Assay Params) – Edit 2 Pt Rate Additional Parameters**

Initial Abs. Limits: -0.20 to 3.50  
Second Abs. Limits: -0.20 to 3.50  
Antigen Excess Factor: 9.0

**DRI Propoxyphene Assay – Semiquantitative  
Ortho Clinical Diagnostics VITROS® XT 7600 System, VITROS® 5600 System, and  
VITROS® 4600 System Parameters, *continued***

**Edit Protocol Parameters**

Step	Volume	Pack ID	Seconds	Wavelength
1. Reagent	100 µL	UDxx /A		
2. Incubation			0.0	
3. Sample	8			
4. Incubation			304.0	
5. Reagent	100 µL	UDxx /B		
6. Incubation			38.0	
7. Read				340 nm
8. Incubation			76.0	
9. Read				340 nm

**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>0</u>
2	<u>1.0</u>	<u>0.20</u>	Cal Value: <u>150</u>
3	<u>1.0</u>	<u>0.20</u>	Cal Value: <u>300</u>
4	<u>1.0</u>	<u>0.20</u>	Cal Value: <u>500</u>
5	<u>1.0</u>	<u>0.20</u>	Cal Value: <u>1000</u>

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

Monotonicity: Increase

Max Resp High: 3.00

Min. Resp. High: 3.00

Cal Fit Goodness Limit: 0.990

Max Resp. Low: -3.00

Min Resp. Low: -3.00

**Edit Triple Read Parameters**

	Reportable Conc.	Triple Read Limit
Reportable Min.:	<u>0</u>	<u>24.00</u>
Critical Conc.:	<u>300</u>	<u>8.0</u> %
Reportable Max.:	<u>1000</u>	<u>8.0</u> %

**Comments: N/A**

---

© 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*