



# Liquichek™ Qualitative Urine Toxicology Control



## Liquichek™ Qualitative Urine Toxicology Control

A liquid control designed for use with qualitative rapid urine drug screen panels, including point-of-care Drugs-of-Abuse test kits like Bio-Rad TOX/See™, Biosite Triage® and other popular test kits. The package insert includes qualitative results (+/-) for popular kits and recovery values by gas chromatography (GC) and other quantitative methods.

- Negative and positive levels
- Human urine based
- 3 year shelf life at 2–8°C
- 30 day open-vial stability at 2–25°C



### Analyte Target Values

Analyte	Negative (ng/mL)	Positive (ng/mL)
<b>Catalog Number</b>	454	455
<b>Amphetamines</b>		
d-Amphetamine	<125	2,000
d-Methamphetamine	<125	2,000
3, 4-Methylenedioxymethamphetamine (MDMA)	<125	750
<b>Barbiturates</b>		
Secobarbital	<80	1,000
<b>Benzodiazepines</b>		
Nordiazepam	<50	3,000
Oxazepam	<150	1,000
<b>Cannabinoids</b>		
11-nor-Δ <sup>9</sup> -THC-9-COOH	<3.75	150
<b>Cocaine</b>		
Benzoylcegonine	<30	4,000
Ethanol (mg/dL)	<20	250
Lysergic Acid Diethylamide (LSD)	<0.10	2.0
Methadone	<100	750
Methaqualone	<100	750
<b>Opiates</b>		
Morphine (Free)	<300	3,000
Oxycodone	<60	150
Phencyclidine (PCP)	<5.0	250
Propoxyphene	<150	750
<b>Tricyclic Antidepressants</b>		
Nortriptyline	<10	2,000

Refer to the package insert for performance and stability claims

### Ordering Information

Cat #	Liquichek™ Qualitative Urine Toxicology Control	Quantity
454	Negative Level . . . . .	6 x 3 mL
455	Positive Level . . . . .	6 x 3 mL
454X	Bilevel MiniPak (1 of each level) . . . . .	2 x 3 mL

Additional urine toxicology controls are available.



**Bio-Rad  
Laboratories**

*For further information, please contact the Bio-Rad office nearest you or visit our website at [www.bio-rad.com/toxcontrols](http://www.bio-rad.com/toxcontrols)*

*Clinical  
Diagnostics Group*

**Website** [www.bio-rad.com/qualitycontrol](http://www.bio-rad.com/qualitycontrol) **U.S.** 1-800-2BIO-RAD **Australia** 61-2-9914-2800 **Austria** 43-1-877-8901 **Belgium** 32-9-385-5511 **Brazil** 5521-3237-9400  
**Canada** 1-514-334-4372 **China** 86-21-64260808 **Czech Republic** 420-241-430-532 **Denmark** +45-4452-1000 **Finland** 358-9-804-22-00 **France** 33-1-47-95-60-00  
**Germany** +49-(0)89-318-840 **Greece** 30-210-7774396 **Hong Kong** 852-2789-3300 **Hungary** +36-1-459-6100 **India** 1-800-180-1224 **Israel** 972-3-9636050  
**Italy** +39-02-216091 **Japan** 81-3-6361-7070 **Korea** 82-2-3473-4460 **Mexico** +52 (55) 5488-7670 **The Netherlands** +31-318-540666 **New Zealand** 64-9-415-2280  
**Norway** 47-23-38-41-30 **Poland** 48-22-3319999 **Portugal** 351-21-472-7700 **Russia** 7-495-721-14-04 **Singapore** 65-6415-3188 **South Africa** 27-11-442-85-08  
**Spain** 34-91-590-5200 **Sweden** 46-8-555-127-00 **Switzerland** 41-61-717-95-55 **Thailand** 662-651-8311 **United Kingdom** +44-(0)20-8328-2000



Printed on recycled paper with soy-based inks

BioSite Triage is a registered trademark of BioSite Inc.

© 2010 Bio-Rad Laboratories, Inc. Printed in the USA 04/10 QSD10-023 Q-1613