



Presenting the TOX/See™ Rapid Urine Drug Screen Tests



When every minute counts, TOX/See™ is the answer.





TOX/See™ Rapid Urine Drug Screen Tests

Fast, flexible and convenient drug testing is more important than ever.

	TOX/See™ Results in less than 5 minutes	< vs >		Leading competitor Results in 15 minutes
---	---	--------	---	--

1 Walkaway Flexibility

Test results are stable and readable for a minimum of one hour to accommodate busy laboratory schedules

2 Comprehensive Test Panels

Simultaneous detection of as many as 12 drugs per cassette

3 Rapid Results

Test results are available in less than 5 minutes

4 One-Step Test Protocol

One-step test protocol allows for easy training

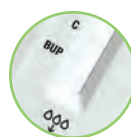
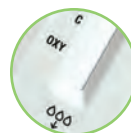
5 Differentiates between Amphetamine, Methamphetamine and Ecstasy in One Step

Provides more specific information to the clinicians for better patient management and saves time and money for the laboratories

6 Specific Tests for Emerging Drugs of Abuse

Oxycodone – From 1994 to 2002, Oxycodone related cases in the US emergency departments increased by 450%. About 70% of these cases involved abuse of Oxycodone along with other drugs*. As this trend continues, the demand for prescription monitoring will multiply.

Buprenorphine – Buprenorphine was recently approved as a substitution treatment for heroin addiction in the USA, following Europe, Australia and Singapore**. As the number of buprenorphine users increases, it will become more important to monitor treatment compliance and efficacy.



* The DAWN Report, June 2004

** <http://www.iatdmct.org/buprenorphine.htm>, Tackling Subutex abuse in Singapore
Lee C E, Singapore Med J 2006; 47(11) : 919

TOX/See™ Rapid Urine Drug Screen Tests

Our Tests Feature a Combination of the Following Drugs at These Cutoffs:

Amphetamine	AMP1000 ng/mL	THC	THC50 ng/mL
Opiates	OPI300 ng/mL	Methamphetamine	MET1000 ng/mL
Opiates	OPI20002000 ng/mL	Barbiturates	BAR300 ng/mL
Benzodiazepines	BZO300 ng/mL	Tricyclic Antidepressants	TCA1000 ng/mL
Methadone	MTD300 ng/mL	Ecstasy	MDMA500 ng/mL
Cocaine	COC300 ng/mL	Phencyclidine	PCP25 ng/mL
Oxycodone	OXY100 ng/mL	Propoxyphene	PPX300 ng/mL
Buprenorphine	BUP10 ng/mL			

Ordering Information

Catalog No.	Description	
194-5221	TOX/See™ Drug Screen Test AMP, BAR, BZO, COC, MDMA, MET, MTD, OPI, OXY, PCP, TCA, THC.....	25 tests
194-5225	TOX/See™ Drug Screen Test AMP, BAR, BZO, COC, MDMA, MET, MTD, OPI, PCP, THC.....	25 tests
194-5228	TOX/See™ Drug Screen Test AMP, BAR, BZO, COC, MET, MTD, OPI, PCP, TCA, THC.....	25 tests
194-5207	TOX/See™ Drug Screen Test AMP, BAR, BZO, COC, MDMA, MET, MTD, OPI, TCA, THC.....	25 tests
194-5230	TOX/See™ Drug Screen Test AMP, BAR, BZO, COC, MET, OPI, PCP, TCA, THC.....	25 tests
194-5182	TOX/See™ Drug Screen Test AMP, BAR, BZO, COC, MET, MTD, OPI, THC.....	25 tests
194-5183	TOX/See™ Drug Screen Test AMP, BAR, BZO, COC, MTD, OPI 2000, PCP, THC.....	25 tests
194-5209	TOX/See™ Drug Screen Test AMP, BAR, BZO, COC, MTD, OPI, PCP, THC.....	25 tests
194-5216	TOX/See™ Drug Screen Test AMP, BZO, COC, OPI, THC.....	25 tests
194-5200	TOX/See™ Drug Screen Test AMP, COC, OPI, PCP, THC.....	25 tests
194-5202	TOX/See™ Drug Screen Test AMP, COC, MET, OPI, THC.....	25 tests
194-5233	TOX/See™ Drug Screen Test AMP, MDMA, MET.....	25 tests
194-5217	TOX/See™ Drug Screen Test AMP, MET, OPI.....	25 tests
194-5215	TOX/See™ Drug Screen Test AMP, OPI, THC.....	25 tests
194-5214	TOX/See™ Drug Screen Test AMP, OPI.....	25 tests
194-5184	TOX/See™ Drug Screen Test AMP.....	25 tests
194-5185	TOX/See™ Drug Screen Test BAR.....	25 tests
194-5205	TOX/See™ Drug Screen Test BUP.....	25 tests
194-5186	TOX/See™ Drug Screen Test BZO.....	25 tests
194-5188	TOX/See™ Drug Screen Test COC.....	25 tests
194-5224	TOX/See™ Drug Screen Test MDMA.....	25 tests
194-5190	TOX/See™ Drug Screen Test MET.....	25 tests
194-5191	TOX/See™ Drug Screen Test MTD.....	25 tests
194-5192	TOX/See™ Drug Screen Test OPI.....	25 tests
194-5203	TOX/See™ Drug Screen Test OXY.....	25 tests
194-5194	TOX/See™ Drug Screen Test PCP.....	25 tests
194-5204	TOX/See™ Drug Screen Test PPX.....	25 tests
194-5234	TOX/See™ Drug Screen Test TCA.....	25 tests
194-5187	TOX/See™ Drug Screen Test THC.....	25 tests



**Bio-Rad
Laboratories**

For further information, please contact the Bio-Rad office nearest you or visit our website at www.bio-rad.com/diagnostics

*Clinical
Diagnostics Group*

Website www.bio-rad.com/diagnostics **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)5200-0520 **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

© 2010 Bio-Rad Laboratories, Inc. Printed in USA 03/10 DG10-0322 C-243