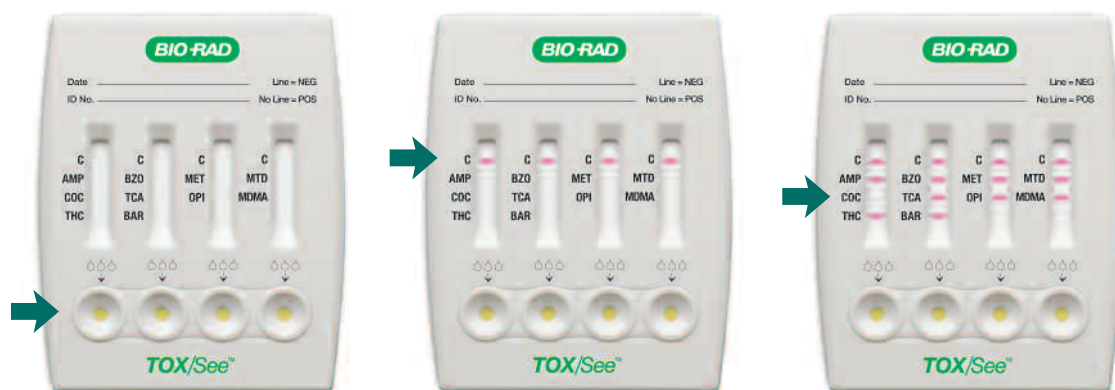


# TOX/See™ Protocol Sheet

## A One-Step Urine Test for Rapid Drug Screen



### 1 Add Urine

- Add 3 drops of urine to each well
- Check control line and read results at 5 minutes after the addition of sample



### 2 Check Control

The presence of a colored line beside the C (control) serves as verification that:

- a sufficient volume of specimen was added
- a proper flow was obtained

### 3 Read Results

- Results remain stable for up to one hour
- Line = negative
- No line = positive

In the example, the test is positive for COC.



## Interpretation



**POSITIVE:** A colored line in the control region (C) and NO COLORED LINE next to the respective drug name



**NEGATIVE:** A colored line in the control region (C) and a COLORED LINE next to the respective drug name. Note: The presence of any line, regardless of intensity, should be interpreted as a negative result



**INVALID:** a control line fails to appear.

- Insufficient specimen or incorrect procedural techniques are most likely reasons for control line failure
- Review procedure and repeat the test with a new test device
- If problem persists, discontinue using the lot immediately and contact Bio-Rad

The example shown is the TOX/See™ 10-drug panel (Catalog number 194-5207).

# TOX/See™ Protocol Sheet

## Intended Use of TOX/See™

- This assay provides only a preliminary analytical test result. A more specific alternate chemical method must be used in order to obtain a confirmed analytical result. The preferred method for confirmation testing and the method established by the Substance Abuse and Mental Health Services Administration (SAMHSA) is gas chromatography/mass spectrometry (GC/MS).
- Clinical consideration and professional judgment should be applied to any drug of abuse test result, particularly when preliminary results are used.
- The use of external positive and negative control material is recommended to test each new lot or shipment of product and every 30 days thereafter.
- This protocol sheet is only for general instructions only. The number of drugs and strips may vary depending on which product has been ordered. For specific information, please refer to the package insert.

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

Amphetamine	AMP	1000 ng/mL	Cocaine	COC	300 ng/mL	Barbiturates	BAR	300 ng/mL
Opiates	OPI	300 ng/mL	Oxycodone	OXY	100 ng/mL	Tricyclic Antidepressants	TCA	1000 ng/mL
Opiates	OPI2000	2000 ng/mL	Buprenorphine	BUP	10 ng/mL	Ecstasy	MDMA	500 ng/mL
Benzodiazepines	BZO	300 ng/mL	THC	THC	50 ng/mL	Phencyclidine	PCP	25 ng/mL
Methadone	MTD	300 ng/mL	Methamphetamine	MET	1000 ng/mL	Propoxyphene	PPX	300 ng/mL

## Ordering Information

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



**Bio-Rad  
Laboratories**

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

*Clinical  
Diagnostics Group*

Website [www.bio-rad.com/diagnostics](http://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 36-1-455-8800 India 91-124-4029300 Israel 972-3-9636050 Italy +39-02-216091 Japan 81-3-6361-6902 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