



Liquichek™ Opiate Control



Liquichek™ Opiate Control

A liquid, third party control containing eight drugs or drug metabolites to monitor the precision of urine toxicology testing procedures. This control has been tested by GC/MS, LC/MS/MS and other methods to confirm the recovery of drug quantities added.

- Human urine
- 3 year shelf-life when stored frozen at -20°C to -70°C
- 30 day open-vial stability at 2–8°C
- Available in two levels, containing drugs or drug metabolites at concentrations above and below most assay cutoffs
- Representative values provided for screening and confirmatory methods



Gravimetric Reference Values

Analyte	Level 1 (ng/mL)	Level 2 (ng/mL)
Catalog Number	416	417
6-Monoacetylmorphine (6-MAM, Heroin Metabolite)	7.5	12.5
Buprenorphine	5	15
EDDP (Methadone Metabolite)	50	150
Fentanyl	0.25	0.75
Meperidine*	100	300
Methadone	150	450
Oxycodone	50	150
Tramadol*	100	300

*International use only. Not for diagnostic use in the United States. Refer to the package insert for performance and stability claims

Ordering Information

Cat #	Liquichek™ Opiate Control	Quantity
416	Level 1	6 x 5 mL
417	Level 2	6 x 5 mL
415X	Bilevel MiniPak	2 x 5 mL

Additional urine toxicology controls are available.



This control is featured in the Unity™ Interlaboratory Program. Participation in the Unity™ Interlaboratory Program can help improve the quality of your laboratory test results and help you meet regulatory and accreditation requirements worldwide. You may also subscribe to one of our Unity™ QC Data Management Solutions designed to improve the effectiveness of your statistical process control. Ask your Bio-Rad Account Representative for more information.

Discover the power of Unity™ Solutions at www.QCNet.com.

BIO-RAD

**Bio-Rad
Laboratories**

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

*Clinical
Diagnostics Group*

Website 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