



Liquichek Opiate Control

BIO-RAD

Liquichek Opiate Control



Targeted for analytes tested in pain management or opioid treatment and compliance programs.

- Contains eight drugs or drug metabolites, at concentrations above and below assay cutoffs
- This control has been tested by GC/MS, LC/MS/MS and other methods to confirm the recovery of drug concentration
- Representative values for screening and confirmatory methods provided
- Human urine
- 3 year shelf-life at -20°C to -70°C
- 30 day open-vial stability at 2-8°C

Gravimetric Reference Values

Analyte (ng/mL)	Level 1	Level 2
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.

EQAS

An independent, external assessment of performance in comparison to your peers.

QCNet.com/eqas

Independent QC

Ongoing, proactive, unbiased daily QC that helps identify errors as they occur or begin to trend.

QCNet.com/independentqc

Unity

QC Data Management tools that help you create a strategy to reduce risk and streamline QC workflow.

QCNet.com/datamanagement

Ordering Information

Catalog No. Description

12005786	Level 1	6 x 5 mL
12005787	Level 2	6 x 5 mL
12002594	Level 1 MiniPak	1 x 5 mL
12002595	Level 2 MiniPak	1 x 5 mL

Refer to myinserts.com or the package insert of currently available lots for specific analyte and stability claims.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

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

Clinical
Diagnostics Group

Website www.bio-rad.com/qualitycontrol U.S. 1-800-224-6723 Australia +61(2)9914-2800 Austria +43-1-877-8901 Belgium +32(3)710-53-00 Brazil +55(31)3689-6600 Canada 1-514-334-4372 China +86 21 6169 8500 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 +1800-180-1224 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 +7-495-721-1404 Singapore +65-6415-3170 South Africa +27-11-442-85-08 Spain +34-91-590-5200 Sweden +46-8-555-127-00 Switzerland +41(0)26-674-55-05/06 Taiwan +886-2-2578-7189 Thailand +662-651-8311 United Kingdom +44(0)20-8328-2000

LIQUICHEK is a trademark of Bio-Rad Laboratories, Inc. in certain jurisdictions.

© 2019 Bio-Rad Laboratories, Inc. Printed in USA QSD19-049 Q-1627 Rev. 05/2019

Printed on recycled paper with soy-based inks