

ReadiUse™ Viral RNA Lysis BufferCatalog number: 60015, 60016
Unit size: 50 mL, 500 ML

Component	Storage	Amount (Cat No. 60015)	Amount (Cat No. 60016)
ReadiUse™ Viral RNA Lysis Buffer	Refrigerated (2-8 °C). Minimize light exposure	50 mL	500 ML

OVERVIEW

Current demands for RT-PCR testing of SARS-CoV-2 have led to a worldwide shortage of RNA extraction kits. A key missing component required by extraction kit manufacturers and hospitals is the RNA Lysis buffer. This ReadiUse™ Viral RNA Lysis Buffer might be used as an alternative for Roche MagNA Pure 96 External Lysis Buffer (Code: 06374913001), Biomérieux NucliSENS Lysis Buffer (Code: 280134) and Qiagen Buffer RNAeasy Lysis Buffer (Code: 79216).

PREPARATION OF WORKING SOLUTION**ReadiUse™ Viral RNA Lysis Buffer working solution**

1. Add 0.5 mL of β -Mercaptoethanol to 50 mL of ReadiUse™ Viral RNA Lysis Buffer.

Note Store the RNA Lysis Buffer containing β -Mercaptoethanol at 4 °C.

SAMPLE EXPERIMENTAL PROTOCOL

The following protocol serves only as a guideline and should be optimized to suit your specific experimental needs. The following protocol was optimized for use with at least 1.5×10^3 cells to a maximum of 5×10^6 cells per purification.

1. Wash the cells once with sterile 1X PBS.
2. Add 100 μ L of ReadiUse™ Viral RNA Lysis Buffer working solution to the washed cells. Mix by pipetting.
3. Purify extracted RNA from the previous step using RNA purification method of choice.

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.