



Viral Inactivation Study of SARS-CoV-2 in Lysis Buffer (LBF)

Methods

Lysis Buffer (LBF) and a PBS control buffer were incubated with SARS-COV-2 virus (150 μ L Lysis Buffer (LBF) or PBS and 200 μ L SARS-COV-2 sample as the RNAadvance Viral XP Instructions for Use indicate. The conditions tested were:

- 20 min Room Temperature
- 60 min Room Temperature
- 20 min heat at 60°C

To reduce Lysis Buffer (LBF) cytotoxic effects, 0.1 mL of each sample was diluted into 0.9 mL PBS. 0.5 mL of each diluted sample was added to a 100 kDa molecular weight (MW) cutoff filter (Amicon) and centrifuged at 10,000 rpm for 5 minutes. Flowthrough was discarded and the remaining sample added and centrifuged. 0.5 mL PBS was added to the sample for buffer exchange, which was followed by another centrifugation. After the 2nd centrifugation, 0.2 mL PBS was added to resuspend the sample and 0.1 mL of the sample was used for TCID₅₀ assay.

In addition to TCID₅₀ assay, 0.1 mL of the virus/buffer mixture was transferred into a single well of a 6-well plate for each condition. Due to cytotoxicity observed, an additional TCID₅₀ assay was performed with the culture media in order to determine whether infectious virus could be detected in these wells.

Results

Controls

The TCID₅₀ for PBS control group is 7.94 E4 after Amicon filtration. This is lower than expected and there was likely some sample loss on the Amicon filters.

Heating (in PBS) at 60°C led to undetectable levels of virus.

Lysis Buffer (LBF)

TCID₅₀ assay

For the LBF buffer, there was toxicity at the 10⁻¹ dilution. There was no evidence of virus induced CPE at 10⁻² dilutions and beyond. This equates to 3.16x 10² TCID₅₀/mL. The heat condition showed the same result.

Buffer	PBS	PBS	PBS	LBF	LBF	LBF
Temperature	RT	RT	60 °C	RT	RT	60 °C
Time (min)	20	60	20	20	60	20
TCID ₅₀ /mL	7.94E4	6.31E4	3.70E1	3.16E2*	3.16E2*	3.16E2*

*maximum titer

TCID₅₀ (culture media)

No virus-induced CPE or cytotoxicity was observed in this assay. This equates to < 1 infectious unit present in the initial 0.1 mL inoculums of LBF-SARS-CoV-2 mixtures (10 TCID₅₀/mL).

Conclusions

The conclusion of this assay indicates a $\geq 7,940$ reduction in viral activity, or $\geq 99.987\%$ effective viral inactivation. These results were consistent at both room temperature and 60°C for 20 min in Lysis buffer (LBF). The Initial TCID₅₀ assay indicated viral inactivation could be confirmed at $\geq 99.602\%$ ($1-(316/7.94 \times 10^4)$) and the concluding culture medial TCID₅₀ assay further improved the sensitivity to $\geq 99.987\%$ ($1-(10/7.94 \times 10^4)$). Confirmation of 100% inactivation is limited by the assay sensitivity due to the cytotoxic properties of Lysis Buffer (LBF).



©2020 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. All other trademarks are property of their respective owners.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at [beckman.com](https://www.beckman.com)

AAG-721206.20