

Rodent Models of Oropouche Virus Infection for Antiviral Screening

Gottshall, S.L.,* Robins, S.L., Drakeley, G., Tamhankar, M., Eskandarinezhad, S., Wasko, A., Molestina, R.E., Rashid, S., and Bradford, R.

American Type Culture Collection, Manassas, Virginia, USA. sgottshall@atcc.org



INTRODUCTION

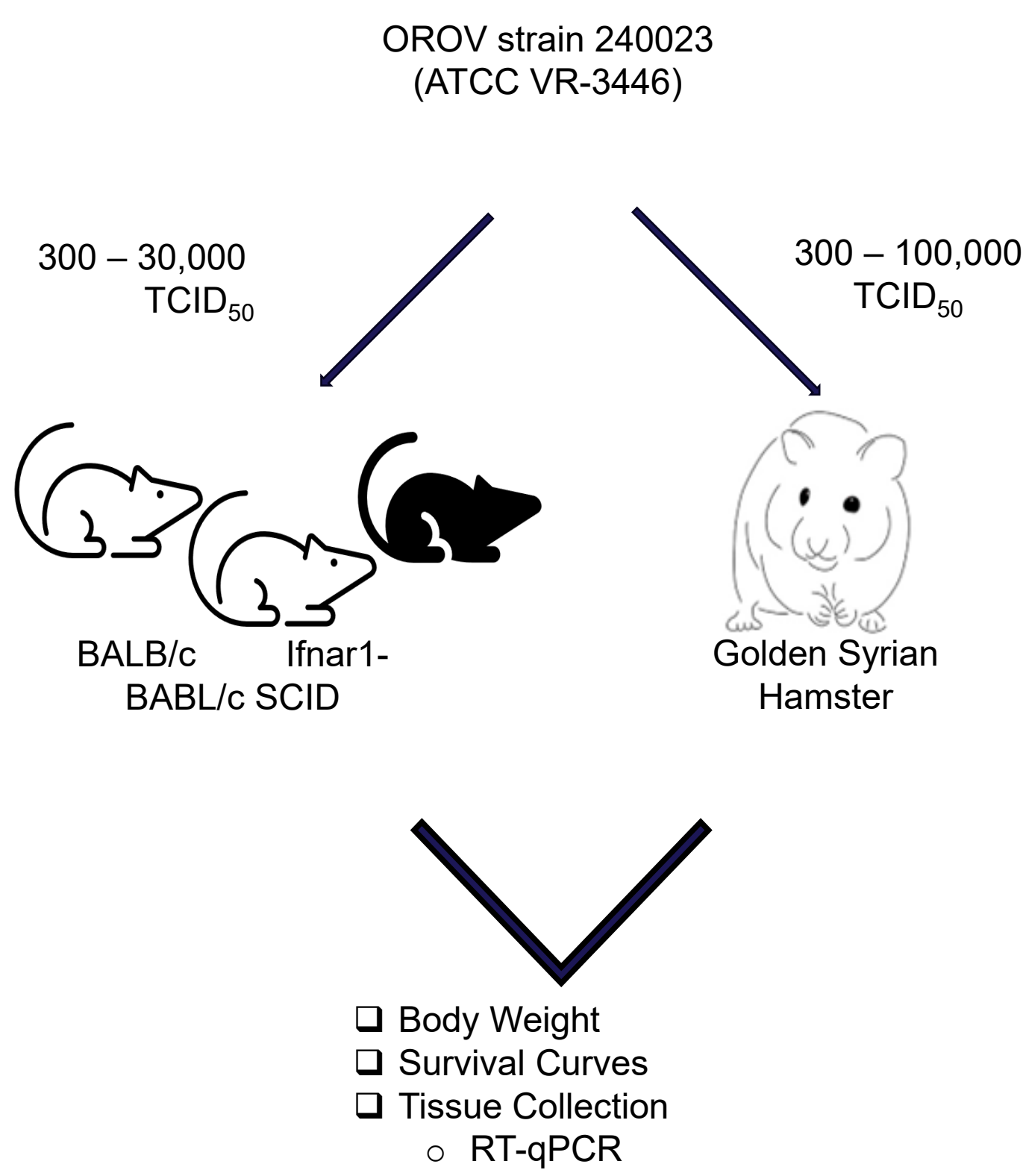
Oropouche virus (OROV) is an emerging public health threat due to its expanding global distribution. The surge in cases is associated with new symptoms, including congenital infection. Genetic analysis revealed this novel symptomology may be associated with emergence of a new natural reassortant. There are no antiviral therapies or vaccines available to treat OROV infection (1-2). Animal models to test the efficacy of antiviral therapies are urgently needed.

EXPERIMENTAL APPROACH

OBJECTIVE: To determine the susceptibility of mice and golden Syrian hamsters to OROV 240023 (ATCC® VR-3446™).

- BALB/c, BALB/c SCID and Ifnar1- mice were infected with 0 - 30,000 TCID₅₀ OROV (N=4/group).
- Golden Syrian hamsters were infected with 0 - 100,000 TCID₅₀ OROV (N=4/group).
- Body weight and clinical observations were performed daily.
- Euthanasia was performed when animals reached predefined terminal endpoints.
- Tissues, including liver, spleen, brain, and serum were collected for viral RNA extraction and RT-qPCR of viral RNA.

Fig. 1: OROV Infection in Mice and Hamsters



RESULTS

Fig. 2. Gross Liver Pathology in OROV Infected Mice



REFERENCES

- Gunter, K., et al. (2024). A reporter Oropouche virus expressing ZsGreen from the M segment enables pathogenesis studies in mice. Journal of virology, 98(9), e0089324. <https://doi.org/10.1128/jvi.00893-24>
- Rodrigues, A. H., et al (2011). Oropouche virus experimental infection in the golden hamster (Mesocricetus auratus). Virus research, 155(1), 35–41. <https://doi.org/10.1016/j.virusres.2010.08.009>

RESULTS

Fig. 3. Susceptibility of Ifnar1- Mice to OROV

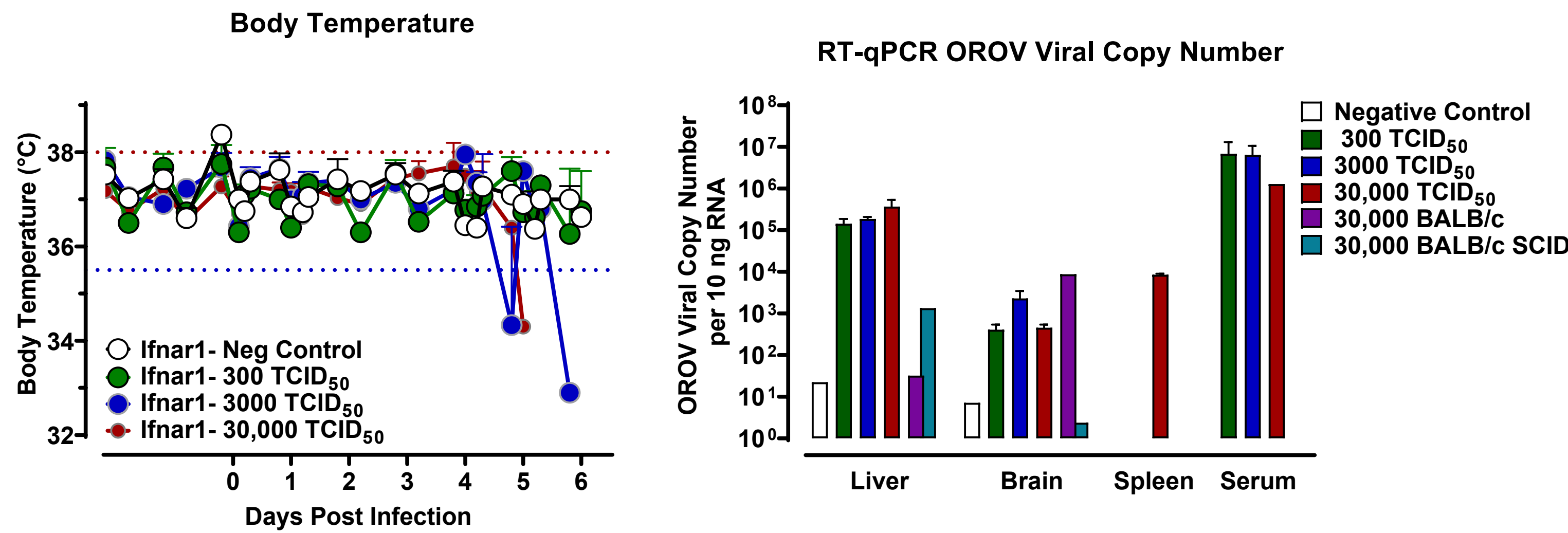
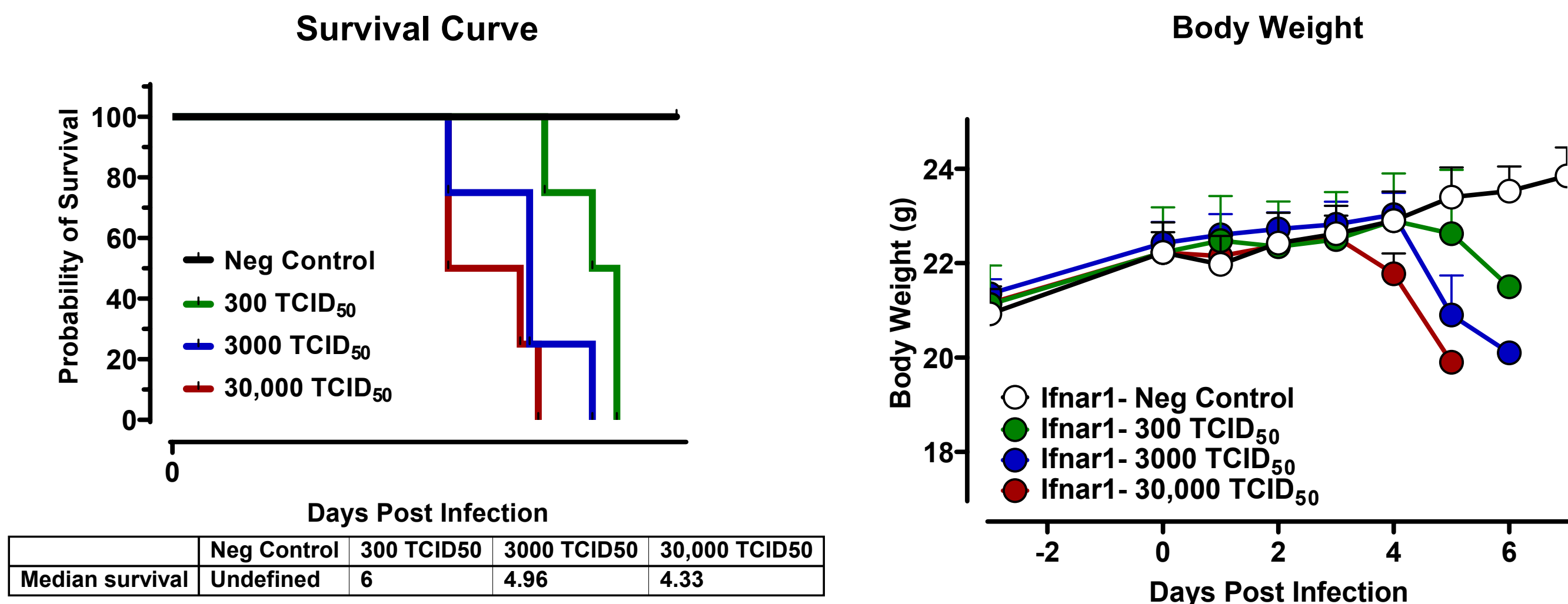
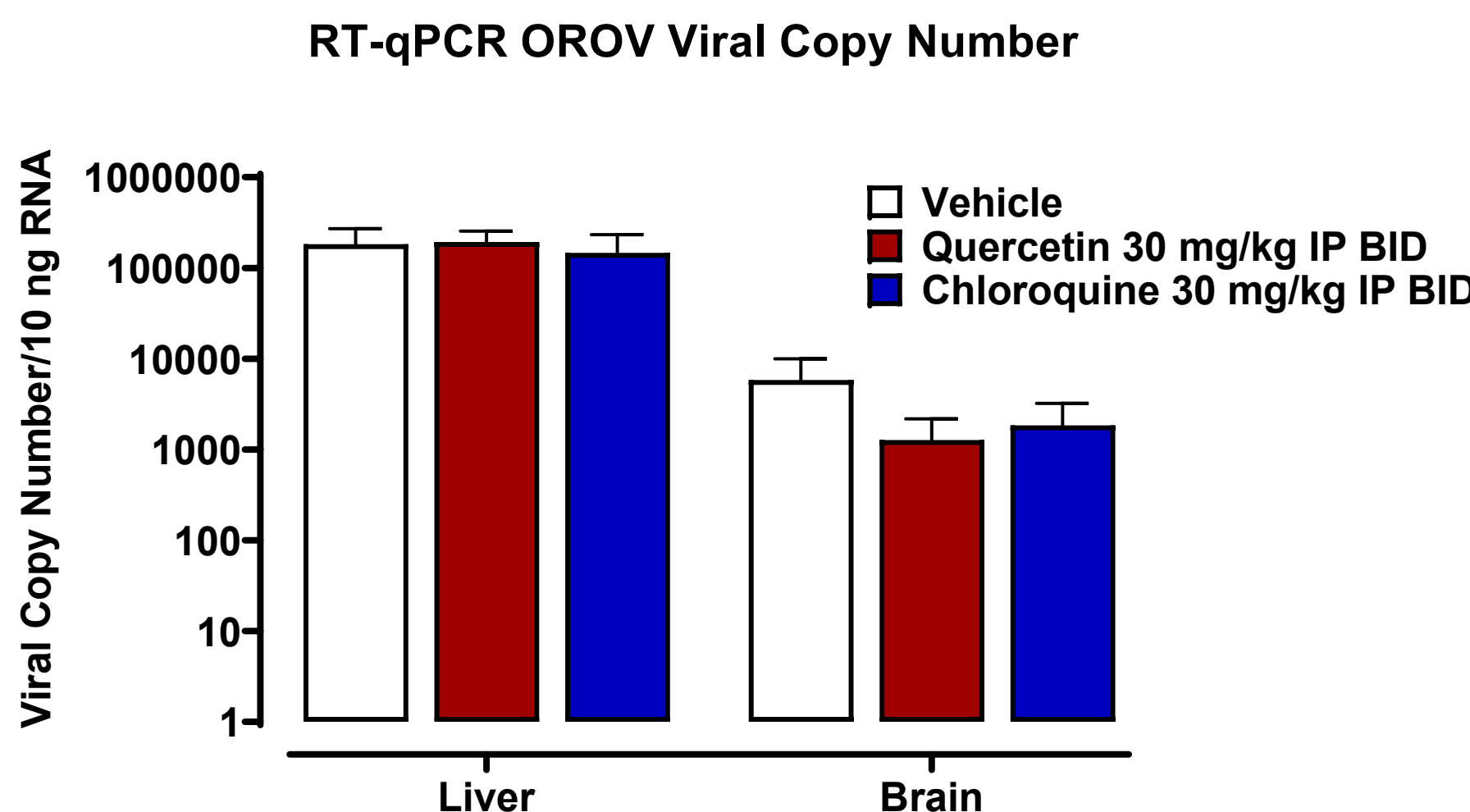
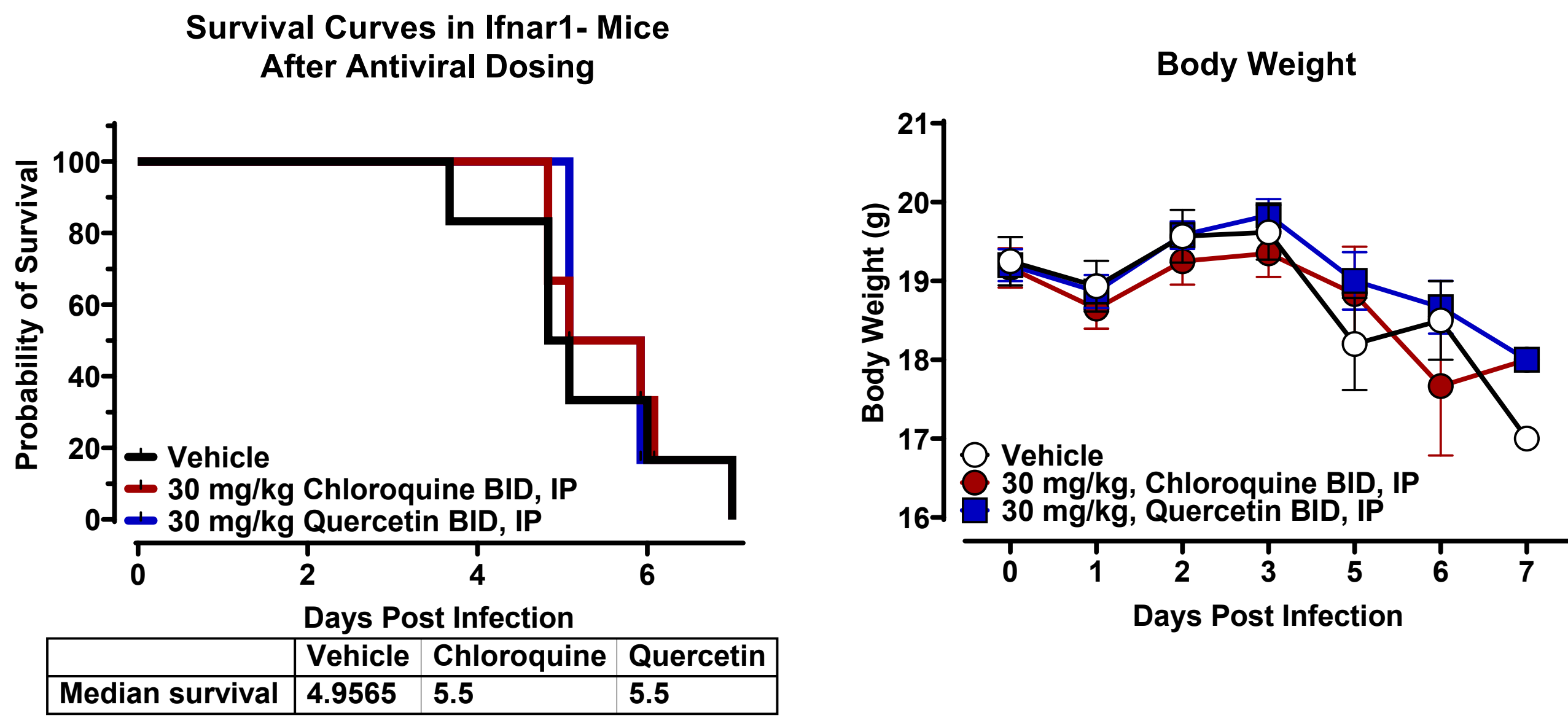


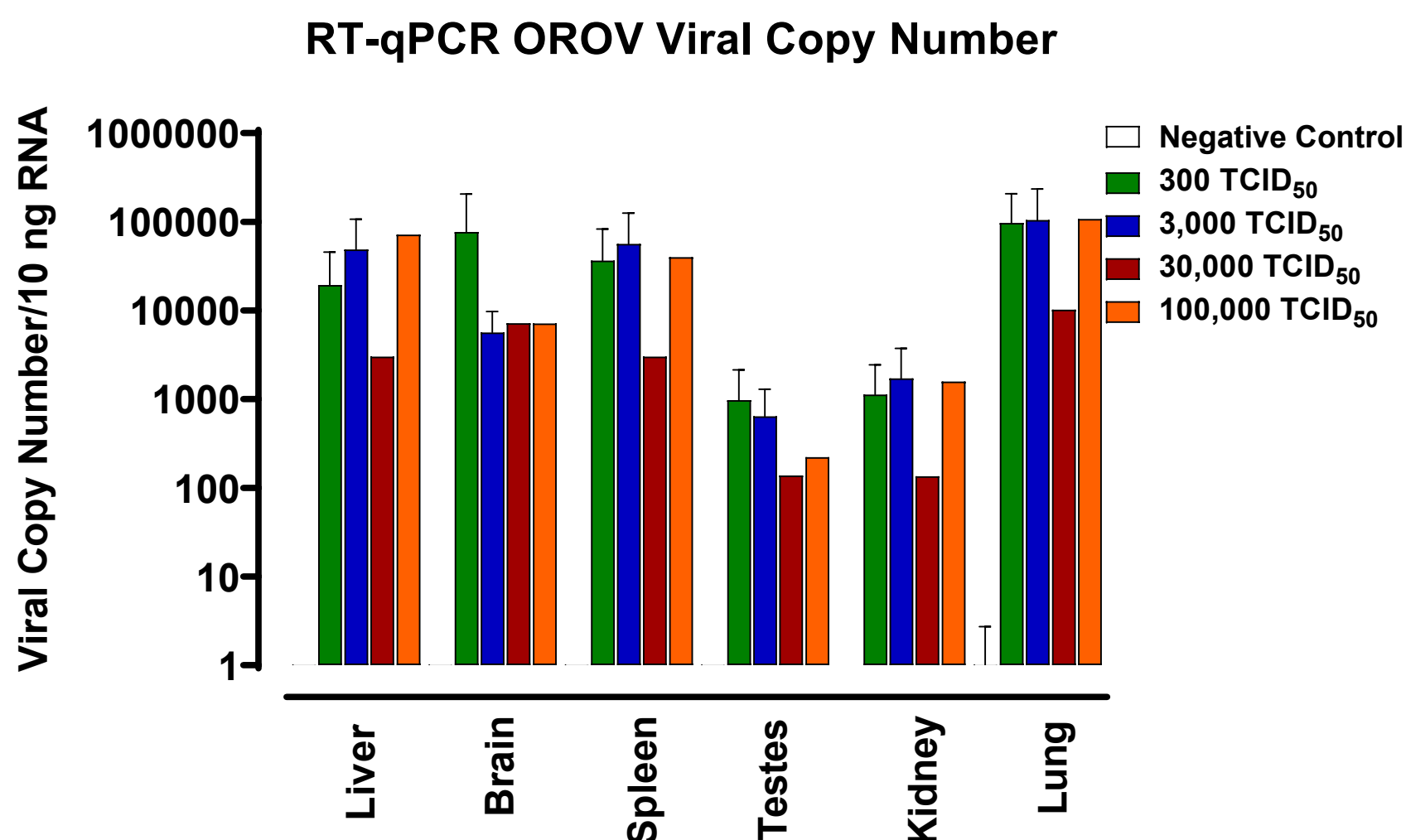
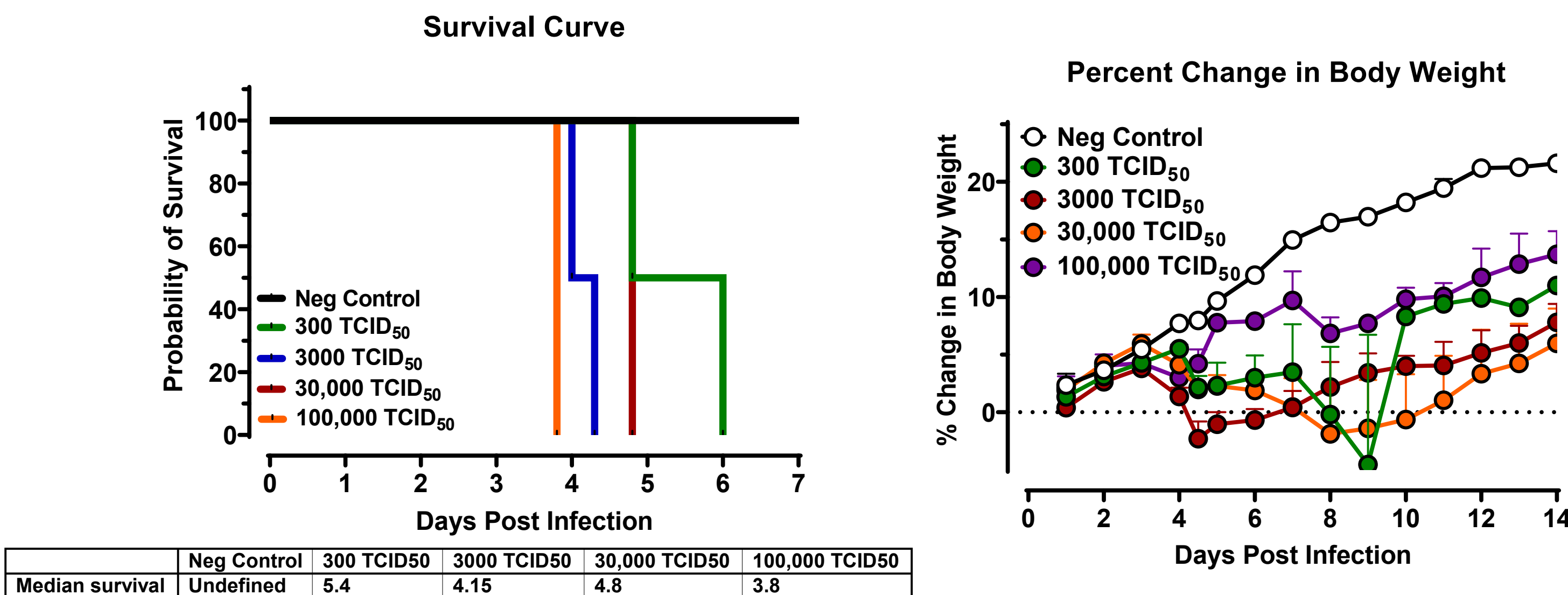
Fig. 4. Antiviral Treatment of OROV Infected Ifnar1- Mice

In a separate study, five-week-old male Ifnar1- mice were infected with 300 TCID₅₀ OROV for 4 h before intraperitoneal dosing of Chloroquine or Quercetin. Antiviral dosing was performed once per day.



RESULTS

Fig. 5. Susceptibility of Golden Syrian Hamsters to OROV



SUMMARY

- OROV infection is 100% lethal in Ifnar1- mice and causes gross liver pathology.
- OROV was rarely lethal in BALB/c (25% at 30,000 TCID₅₀) and BALB/c SCID (25% at 30,000 TCID₅₀) mice.
- Quercetin and Chloroquine showed no antiviral activity against OROV infection in Ifnar1- mice.
- Evidence of OROV infection was detected in 87% of hamsters examined. Lethality was observed in 50% of infected animals. Six hamsters with nonlethal infection showed substantial transient weight loss.
- Of all tissue samples tested, the highest levels of viral RNA were found in the serum of infected mice. Measurable levels of viral RNA were detected in all tissues examined.
- The Ifnar1- mouse model can thus serve as a predictive model in evaluating potential therapies for this emerging disease.

ACKNOWLEDGEMENTS

OROV strain 240023 is available from ATCC® as VR-3446™. This work was performed with funds from the ATCC® Internal Research and Development Program. © ATCC 2019. The ATCC trademark, trade name, any and all ATCC catalog numbers listed in this presentation are trademarks of the American Type Culture Collection unless indicated otherwise.