

Role of the CD47-SIRPa axis in the clearance of *Babesia microti* in vivo



Shamim Mohammad, Robert Molestina*, Ioana Brasov, Susan Gottshall, Sarah Robins, Sujatha Rashid, and Rebecca Bradford.

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

BACKGROUND

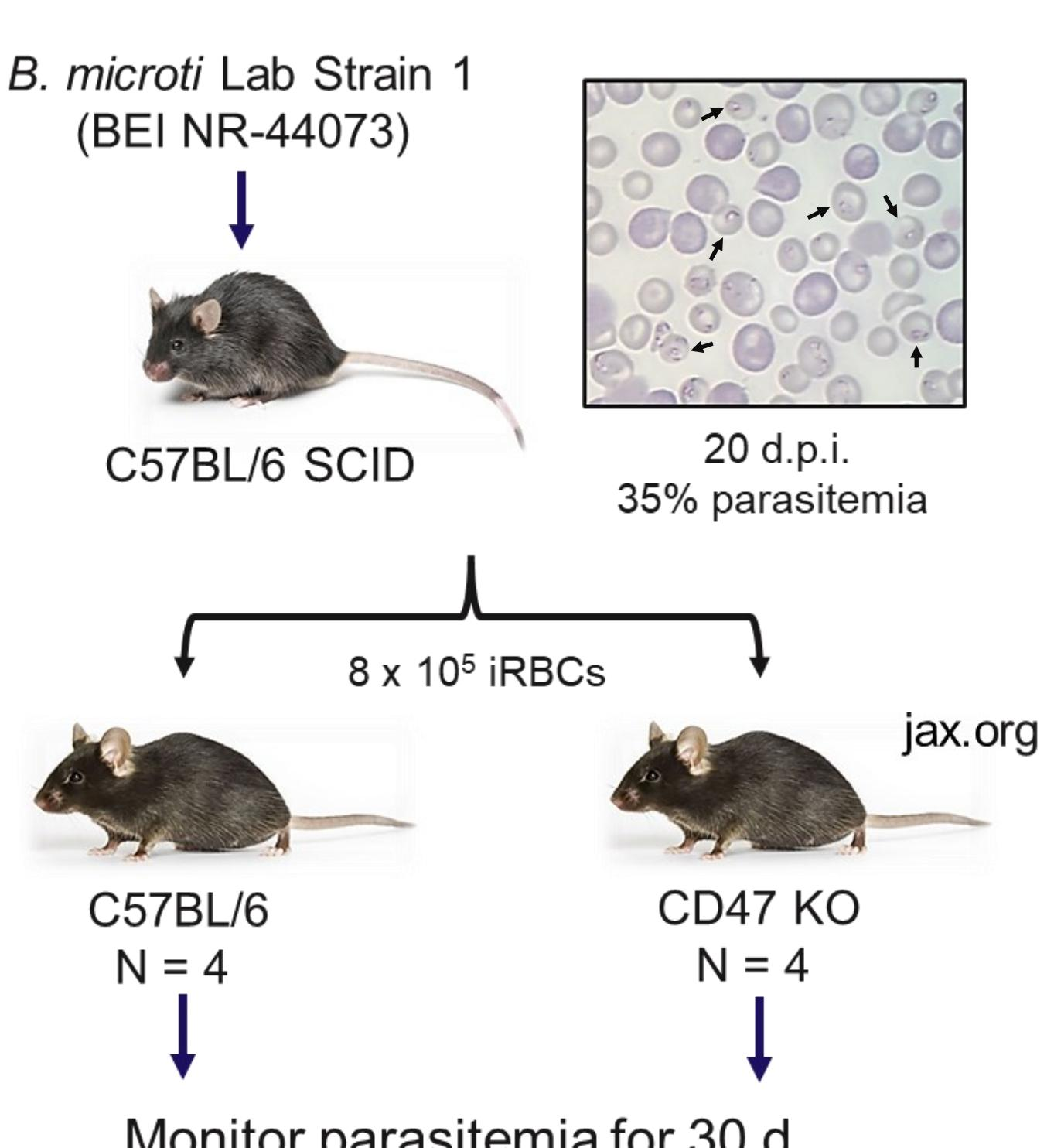
Human babesiosis is an emerging tickborne disease in the US caused by the RBC infecting parasite *Babesia microti*. Infection is primarily transmitted by *Ixodes* ticks, and less frequently by pregnancy or transfusion. Infections are usually asymptomatic or mild among healthy individuals but potentially severe in immunocompromised patients (1). The mechanisms underlying pathogenesis and immune response remain unresolved. In addition, novel therapeutic approaches are needed given the risk for relapsing babesiosis with immunodeficiency (2, 3).

CD47 is expressed on hematopoietic and non-hematopoietic cells and interacts with macrophage signal regulatory protein alpha (SIRPa) and thrombospondin-1 (TSP-1) (4). This interaction induces an antiphagocytic "don't eat me" signal (5, 6).

Hypothesis. *B. microti* preferentially infects young RBCs with high expression of CD47 resulting in protection from phagocytosis. A disruption of CD47 signaling is likely to increase the clearance of *B. microti*-infected RBCs.

EXPERIMENTAL APPROACH

Fig. 3. Infection of CD47 KO Mice

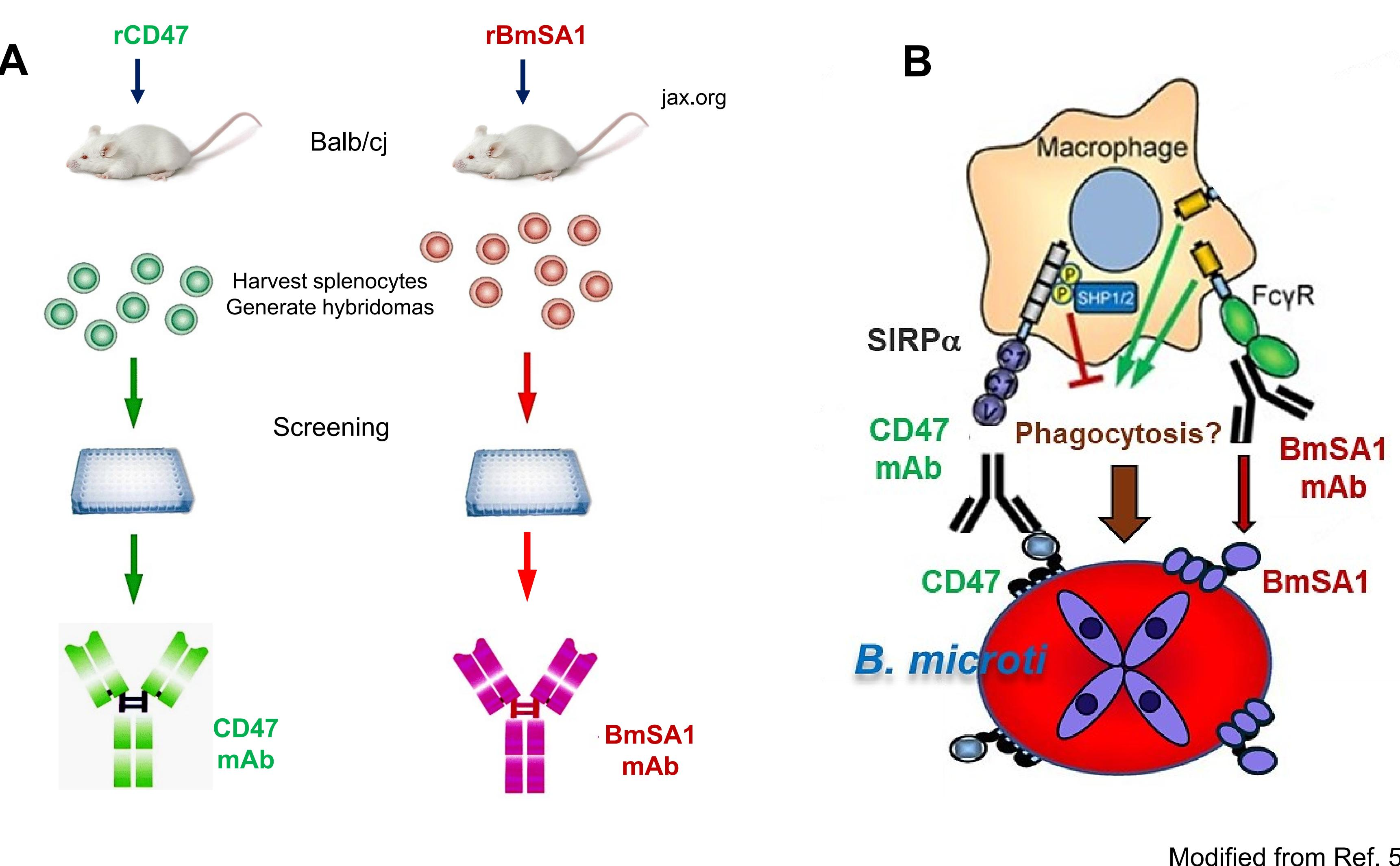


Objective. To investigate the role of CD47 on the growth of *Babesia microti* in mice.

- *B. microti* Lab Strain 1 (BEI NR-44073) was propagated in C57BL/6 SCID mice.
- Eight-week-old female C57BL/6 and CD47 KO (B6.129S7-Cd47^{tm1Fpl}) mice were injected i.p. with 8×10^5 infected red blood cells (iRBCs).
- Blood samples were collected from the tail veins and parasitemia was monitored over a 30-d period of infection by microscopic examination.

- A. In separate experiments, monoclonal antibodies (mAbs) to recombinant CD47 and *B. microti* surface antigen BmSA1 were generated in Balb/c mice.
- B. The utility of these mAbs to block CD47 signaling and promote parasite clearance in susceptible mouse strains is the subject of current research in our laboratory.

Fig. 4. Generation of mAbs to study the role of CD47 in babesiosis



RESULTS

Fig. 5. Susceptibility of CD47 KO mice to *B. microti*

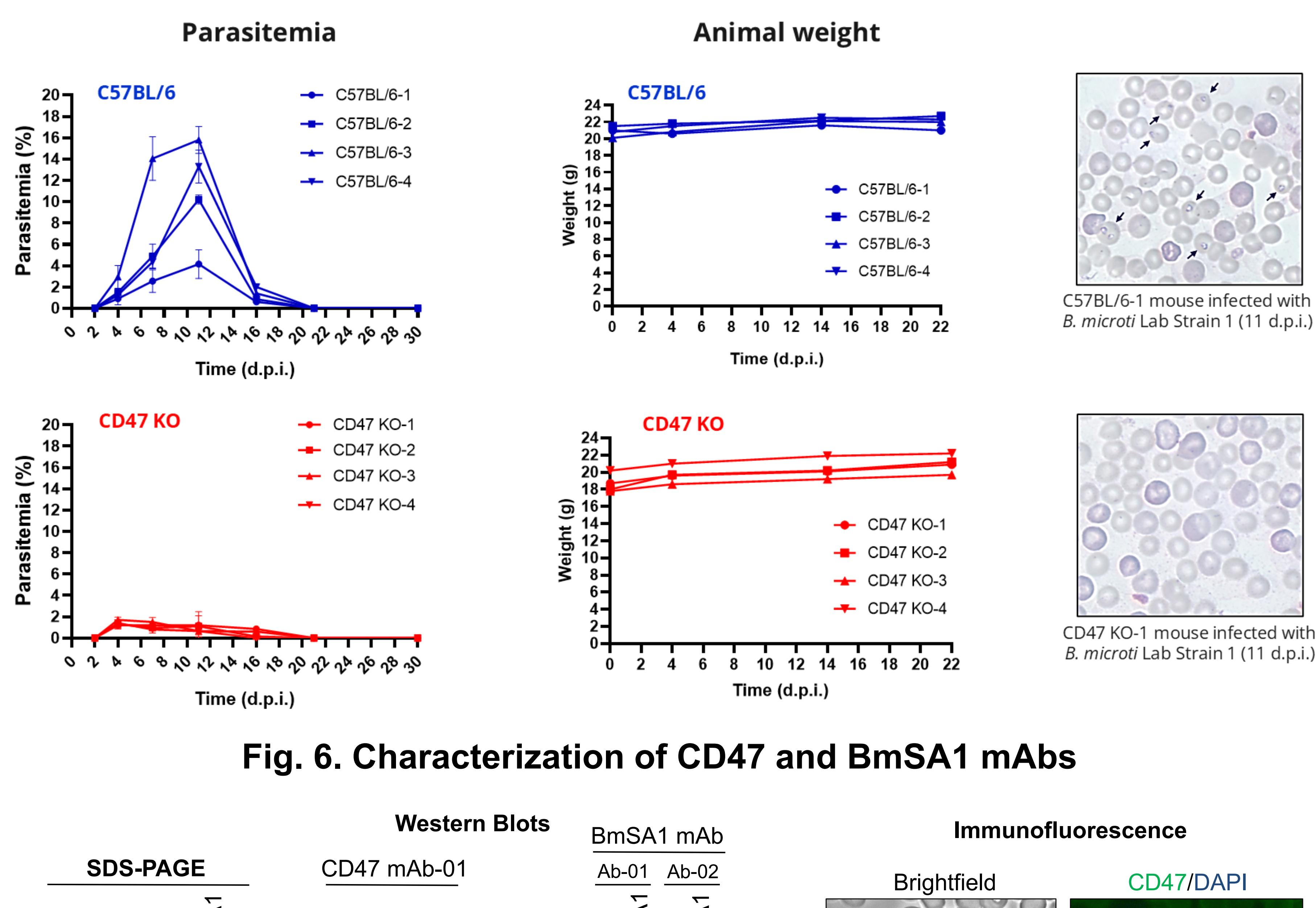
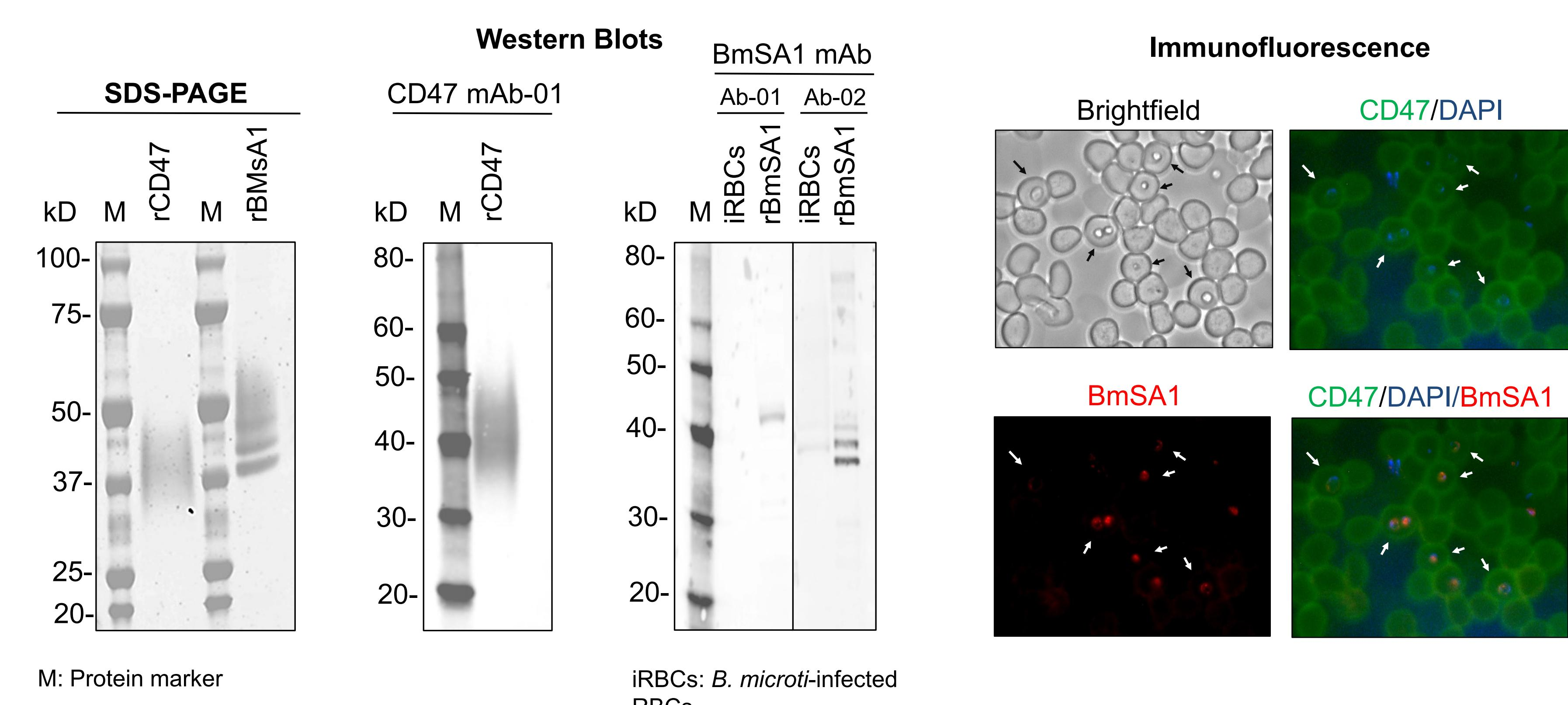


Fig. 6. Characterization of CD47 and BmSA1 mAbs



SUMMARY

- The present study determined the effect of CD47 loss on *B. microti* growth by comparing parasitemia in wild-type C57BL/6 and CD47 KO mice.
- On average, C57BL/6 mice developed 6% parasitemia by day 7 and reached peak parasitemia of 11% by day 11, with infection self-resolving by day 21.
- In contrast, CD47 KO mice developed markedly low parasitemia that on average did not reach >1.5% throughout infection and resolved by day 16.
- We generated and characterized mAbs to CD47 and the *B. microti* antigen BmSA1 for use in subsequent studies in infected SCID mice.
- Future studies will examine the correlation between CD47 level on RBCs and parasite burden, the capacity of *B. microti* to invade CD47 KO RBCs, and the immunological basis underlying the observed CD47-mediated resistance in babesiosis.

REFERENCES

1. Krause, P.J. 2019. Int. J. Parasitol. 49:165-174.
2. Kumar, A., O'Bryan, J., Krause, P.J. 2021. Pathogens 10: 1447.
3. Rogers, R., Krause, P.J., Norris, A.M., Ting M. H. 2023. Clin. Infect. Dis. 76: 741-744.
4. Maute, R., Xu, J., Weissman, I.L. 2022. Immunoncol. Technol. 13: 100070.
5. Kaur, S., Cicalese, K.V., Banerjee, R., Roberts, D.D. 2020. Antib. Ther. 3: 179-192.
6. Burger, P., de Korte, D., van den Berg, T.K., van Bruggen, R. 2012. Transfus. Med. Hemother. 39: 348-352.

ACKNOWLEDGMENTS

Babesia microti Lab Strain 1, NR-44073, was obtained through BEI Resources, NIAID, NIH (www.beiresources.org).

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.