

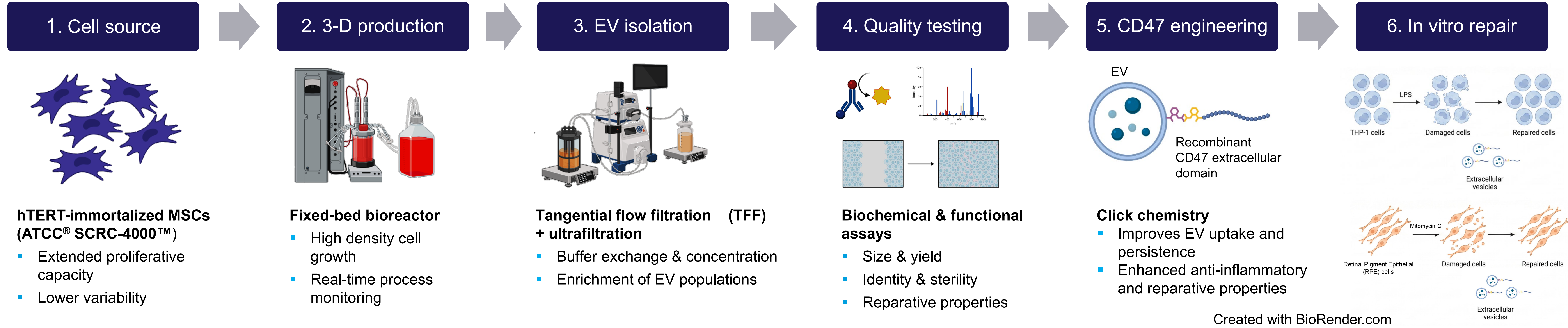
A Scalable Platform for Producing and Surface-Engineering Extracellular Vesicles with Enhanced Reparative Properties

Zachary Cuba, MS;¹ Lenny Godinho, BS;¹ Sujata Choudhury, PhD;¹ David Routenberg, PhD;² Fatah Kashanchi, PhD;³ Heather Couch, PhD;¹ Heather Branscome, PhD^{1,3}
¹ ATCC, Manassas, Virginia, USA, ² Meso Scale Diagnostics, LLC, Rockville, Maryland, USA, ³ George Mason University, Manassas, Virginia, USA

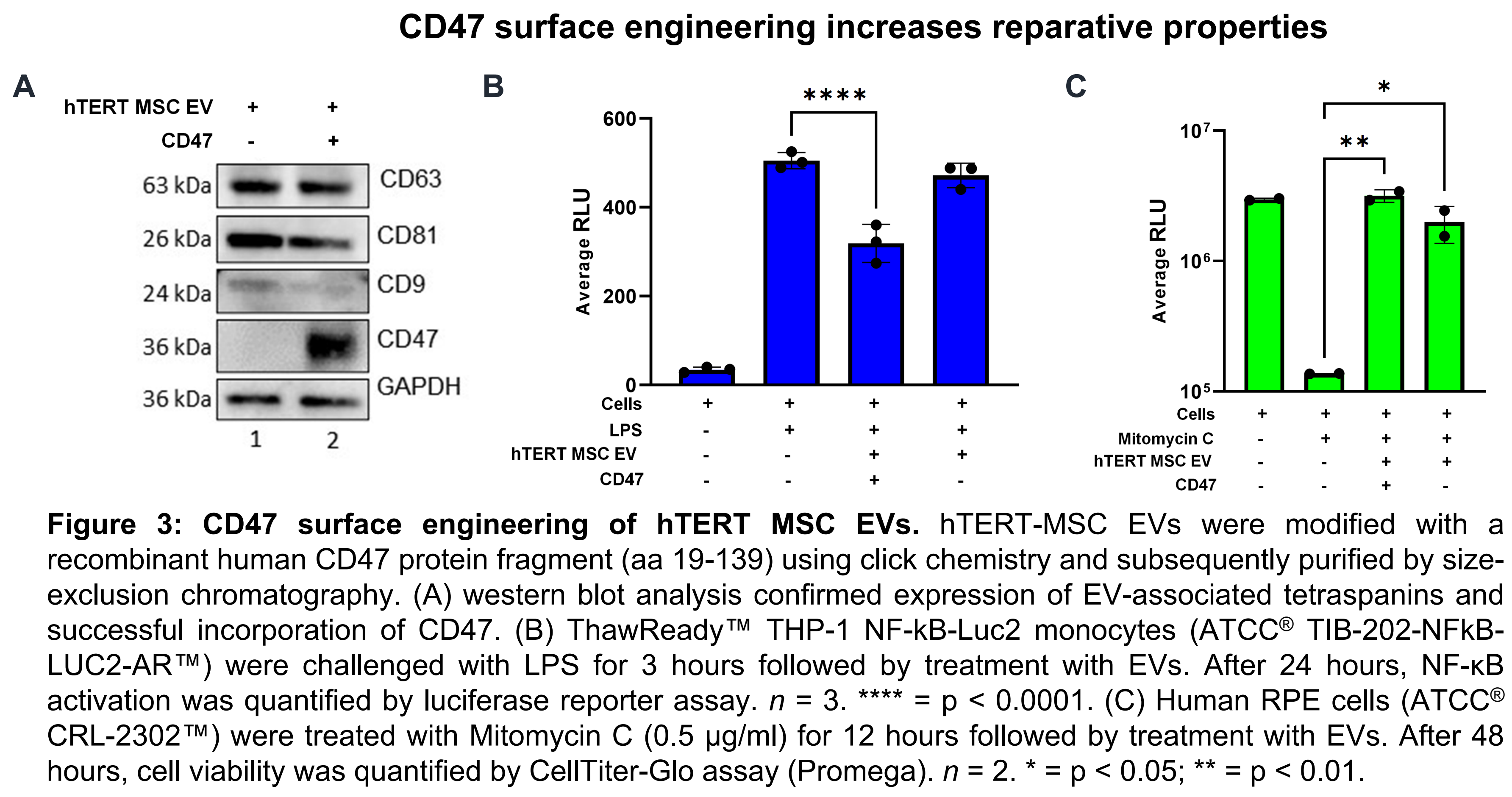
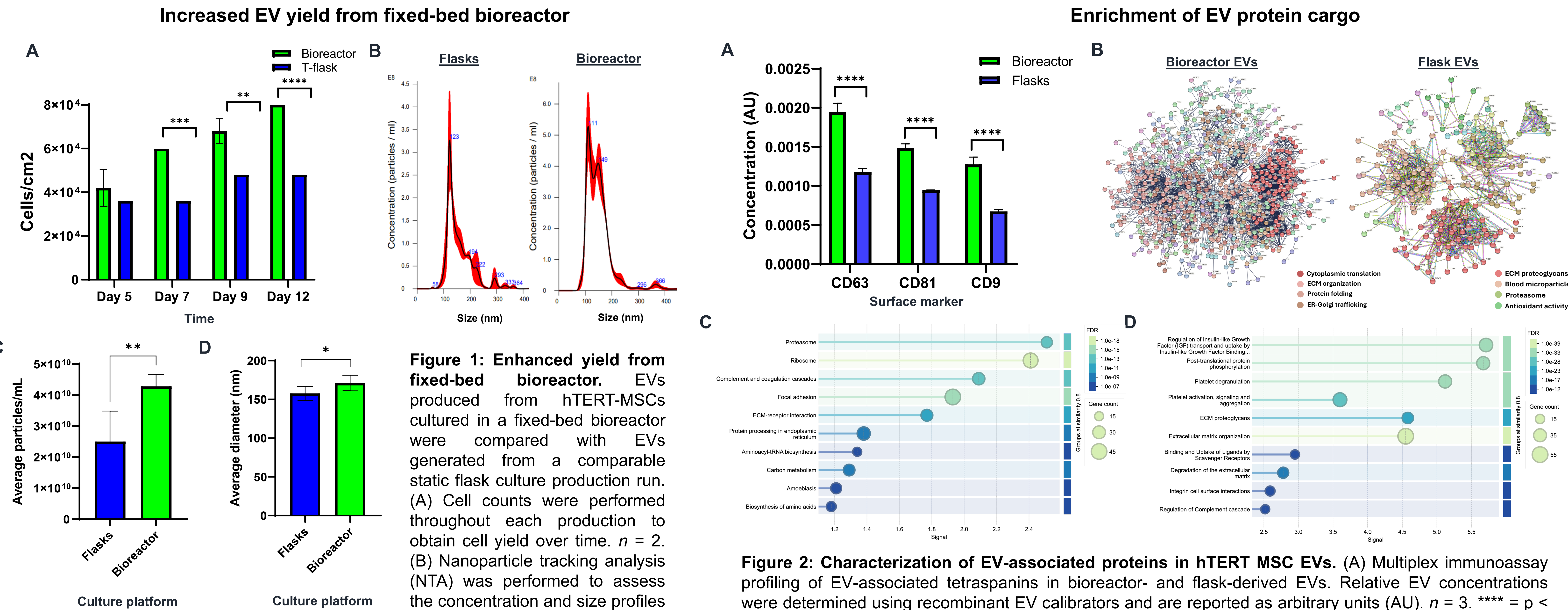
Introduction

- Extracellular vesicles (EVs) are cell-derived nanoparticles that facilitate intercellular communication through the transfer of bioactive cargo.
- Mesenchymal stromal cell (MSC)-derived EVs are widely studied for their immunomodulatory and tissue-supportive biological properties.
- Inflammation and tissue injury are common features of many conditions relevant to military health.
- Broader adoption of EV technologies is limited by challenges in scalable manufacturing, product consistency, and effective delivery to target cells.
- This study combined scalable EV production with CD47 surface engineering, a "don't eat me" signal that may enhance EV persistence and cellular uptake.

Methods



Results



Conclusions

- Scalable production:** Fixed-bed bioreactor culture substantially increased EV yield compared to conventional flask culture.
- Molecular integrity:** Bioreactor-derived EVs retained characteristic EV features and complex protein cargo.
- Biological relevance:** Functional activity was maintained following scale-up and further enhanced through CD47 surface engineering.



Explore
exosomes