

Developing a Microglial HIV-1 Latency Reporter Cells for Investigating Viral Latency and Treatment Strategies



Janice Yoon¹, Helen Navin¹, Scott Anthony², Sameer Tiwari², Yuntao Wu², Sujatha Rashid¹, Rebecca Bradford¹ and Khalid Timani^{1*}

¹ATCC, Manassas, VA 20110 (*Contact: ktimani@atcc.org), ²Center for Infectious Disease Research, George Mason University, Manassas, VA 20110

Abstract

Despite the success of combination antiretroviral therapy (cART) in controlling HIV replication, a definitive cure remains unattainable due to persistent viral reservoirs¹. A major barrier is the establishment of latent HIV within cellular reservoirs, particularly in microglia, the resident macrophages of the central nervous system (CNS)². These cells serve as long-lived viral sanctuaries, contributing to viral persistence, drug resistance, and the pathogenesis of HIV-associated neurocognitive disorders (HAND). Targeting HIV-infected microglia is thus critical for achieving a functional cure.

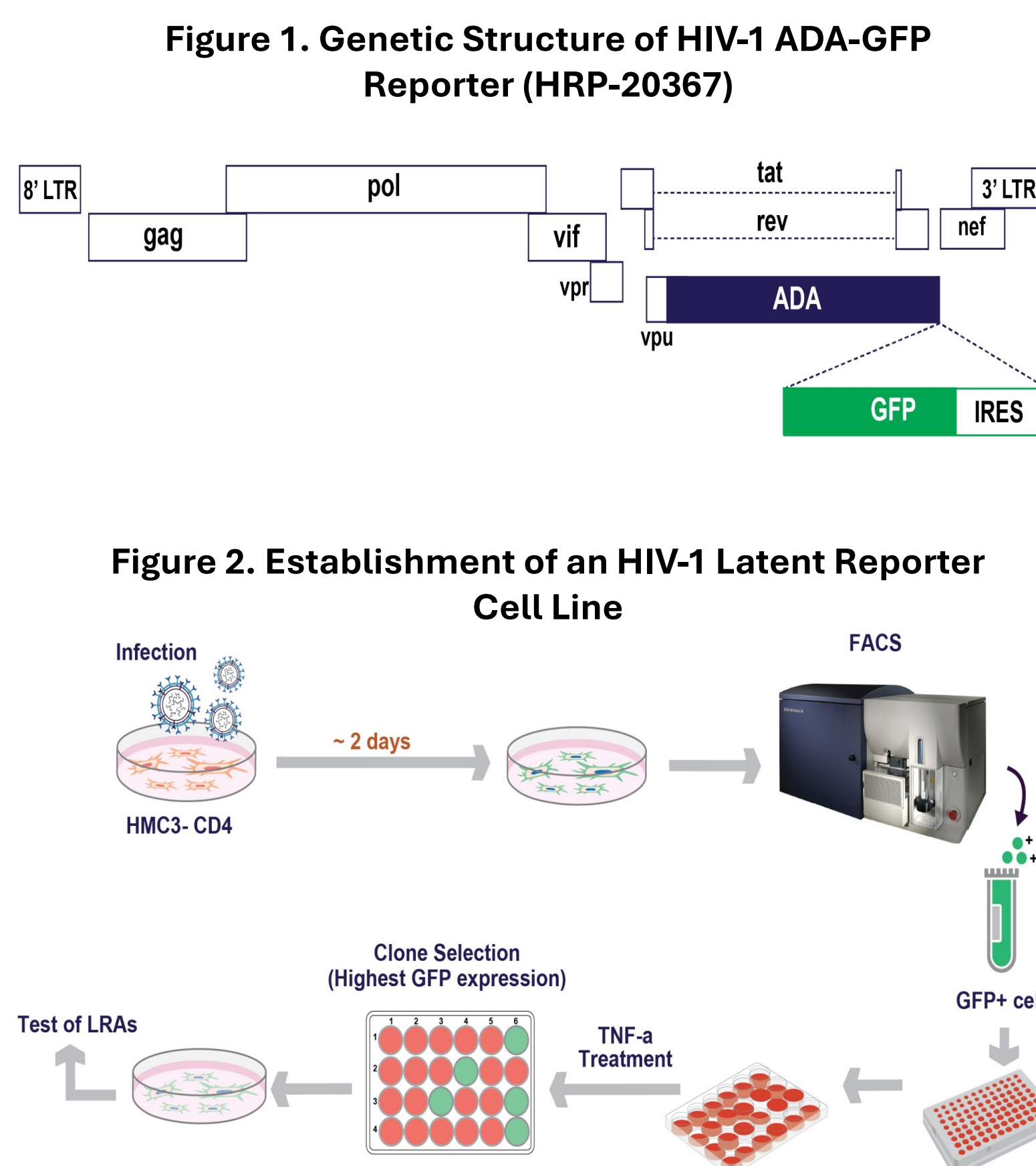
To address this challenge, we aim to develop a human microglial HIV-1 latency reporter cell line using the well-authenticated HMC3 cell line³ (ATCC® CRL-3304). This model will enable mechanistic studies of viral latency in the CNS, facilitate drug screening, and provide a platform to investigate HAND. Furthermore, the latently infected microglia can be incorporated into microglia-containing human brain organoids (MC-HBOs), a 3D in vitro system that recapitulates key features of the human brain, will be a valuable tool to study HIV-CNS interactions, and the efficacy of new therapeutic and curative strategies on the CNS viral reservoir. While we have successfully established the first HIV-1 infected microglia HMC3 cells, current infectivity levels remain low. Ongoing efforts are focused on enhancing infectivity and establishing latent HIV-1, with GFP reporter expression under viral promoter control enabling real-time monitoring of latency reversal and reactivation dynamics.

Methods

NL4.3 ADA-GFP-IRES-Nef virus. An infectious clone (Fig. 1) was used to generate purified virus. Additionally, a dual envelope approach was employed, incorporating VSG to enhance the viral entry. The virus contains the CCR5-tropic ADA Env and co-expresses Nef and GFP from a single bicistronic RNA by IRES.

Overview Establishment HIV-1 latent reporter microglial cell line.

HMC3 stably expressed CD4 cells will be infected with an ADA-GFP reporter. At 2–5 dpi, GFP⁺ cells will be isolated via FACS and plated into 96-well plates. Sorted cells will be monitored for 15–30 days to confirm the establishment of latency, as indicated by diminished GFP expression. Following expansion, The infected cells will be treated with TNF- α to reactivate HIV-1. Clones exhibiting the highest GFP induction (indicating viral reactivation) will be selected, cryopreserved, and subsequently used for screening of latency-reversing agents (LRAs) (Fig.2).

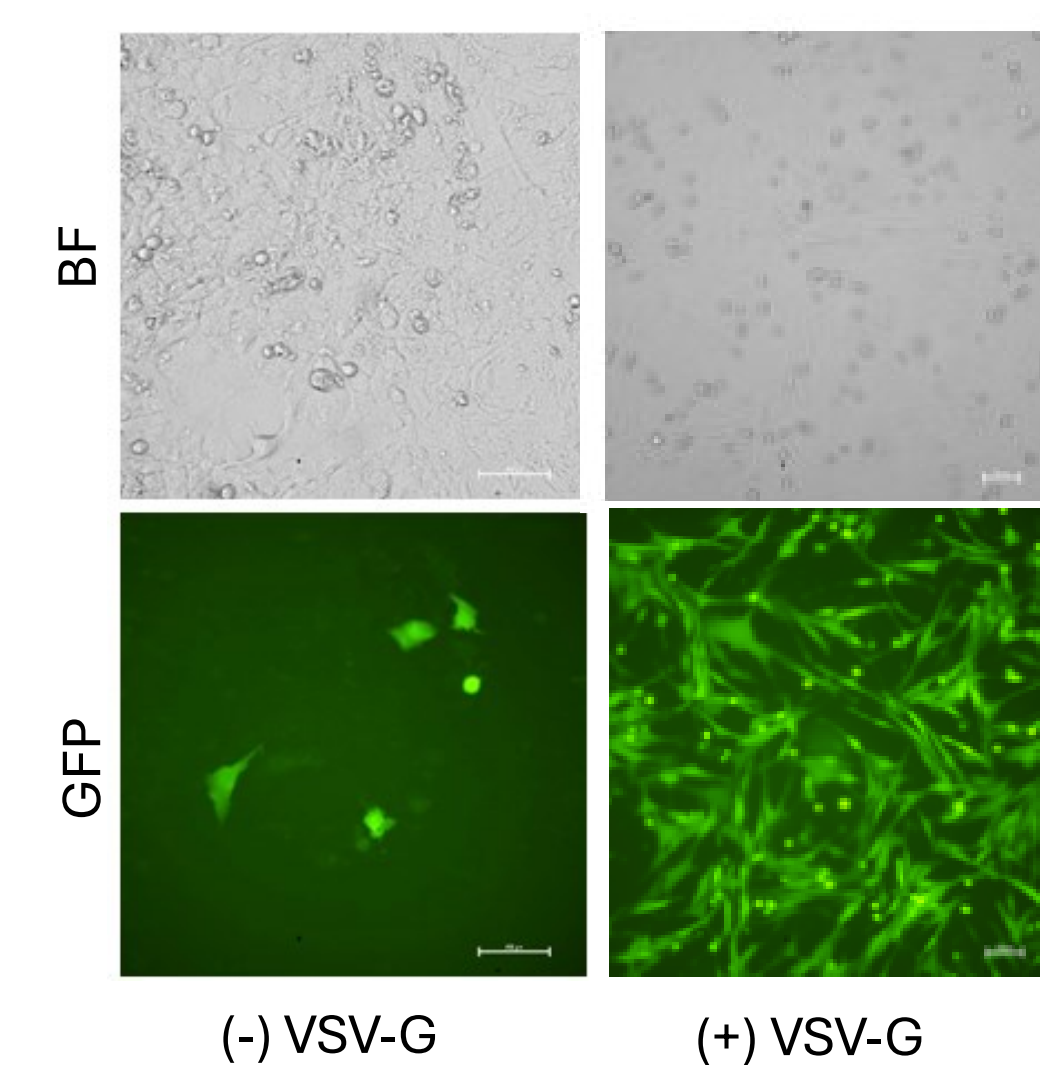


Results

Generation of HMC3-CD4 Stable Cell Line

Due to the lack of the HIV CD4 receptor, HMC3 was engineered for stably expression via lentiviral transduction with a CD4-expressing vector. Several stable clones were generated, and based on moderate CD4 expression levels, a stable cell line (Clone 9) was selected. Flow cytometry analysis confirmed the expression of CD4 as well as the co-receptors CCR5 and CXCR4 in the HMC3-CD4 clone 9 stable cells. Additionally, antibody titration was performed to determine the optimal antibody concentration for cell sorting (Fig 3).

Figure 4. HMC3-CD4 clone 9 infection by HIV-1 ADA-GFP virus



Infection of HIV-1 Reporter Virus in HMC3-CD4 Clone 9

To assess the infectivity of HMC3-CD4 stably expressed cells, the cells were infected with ADA-GFP virus. Interestingly, the cells were successfully infected, as indicated by GFP expression, though the infection level was low (1–2%). To further confirm viral infectivity, we used a dual-envelope virus (+ VSV-G), which showed a significantly higher infection rate (~90%). These data suggest that CD4 expression levels may influence viral infection efficiency (Fig. 4).

Minimal Enhancement of ADA-GFP Infectivity Observed in Higher CD4+ HMC3 subclone

Due to the low infectivity observed in HMC3-CD4 Clone 9 cells, we further sorted these cells based on CD4 expression levels using FACS. The cells were first gated for coreceptors, CCR5⁺/CXCR4⁺, populations and then sorted into six distinct subclones based on CD4 expression (Fig. 5A). Each subclone was infected with ADA-GFP, and infectivity was monitored via GFP expression. Notably, Subclone 9-6, which exhibited the highest CD4 expression, showed only a minimal (~5%) enhancement in infectivity compared to other subclones (Fig. 5B).

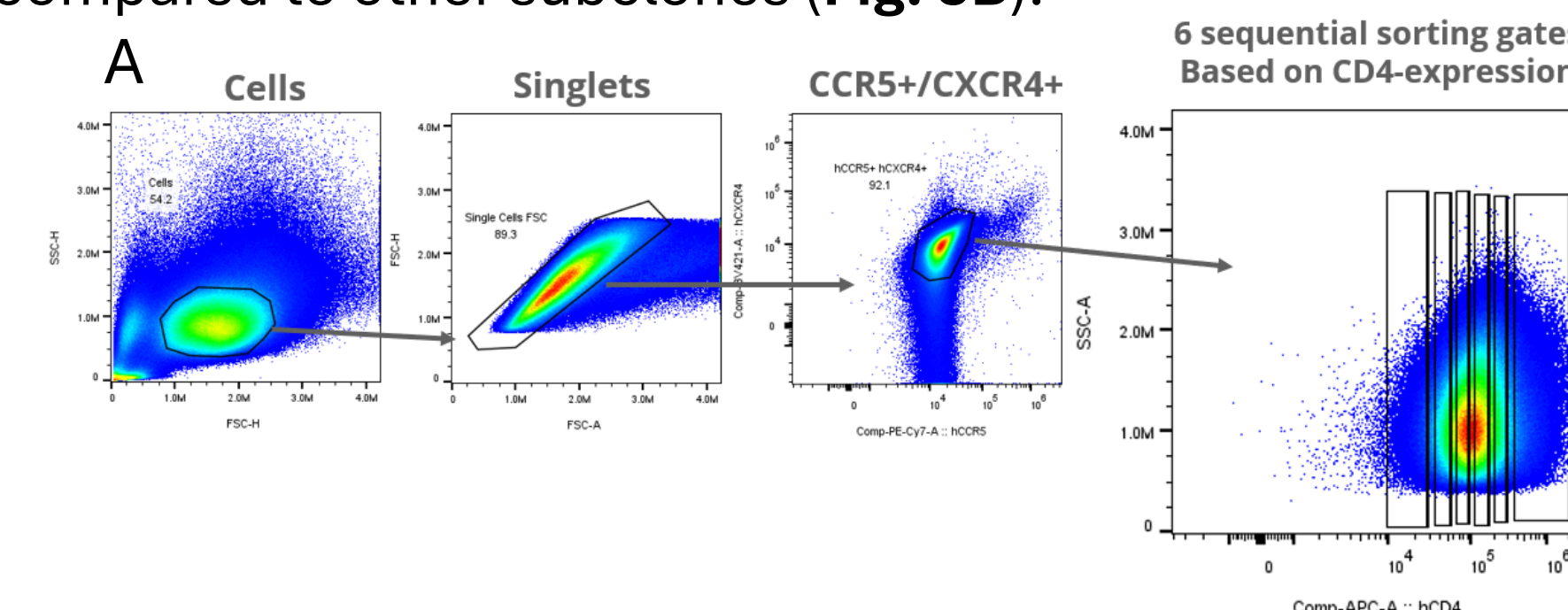
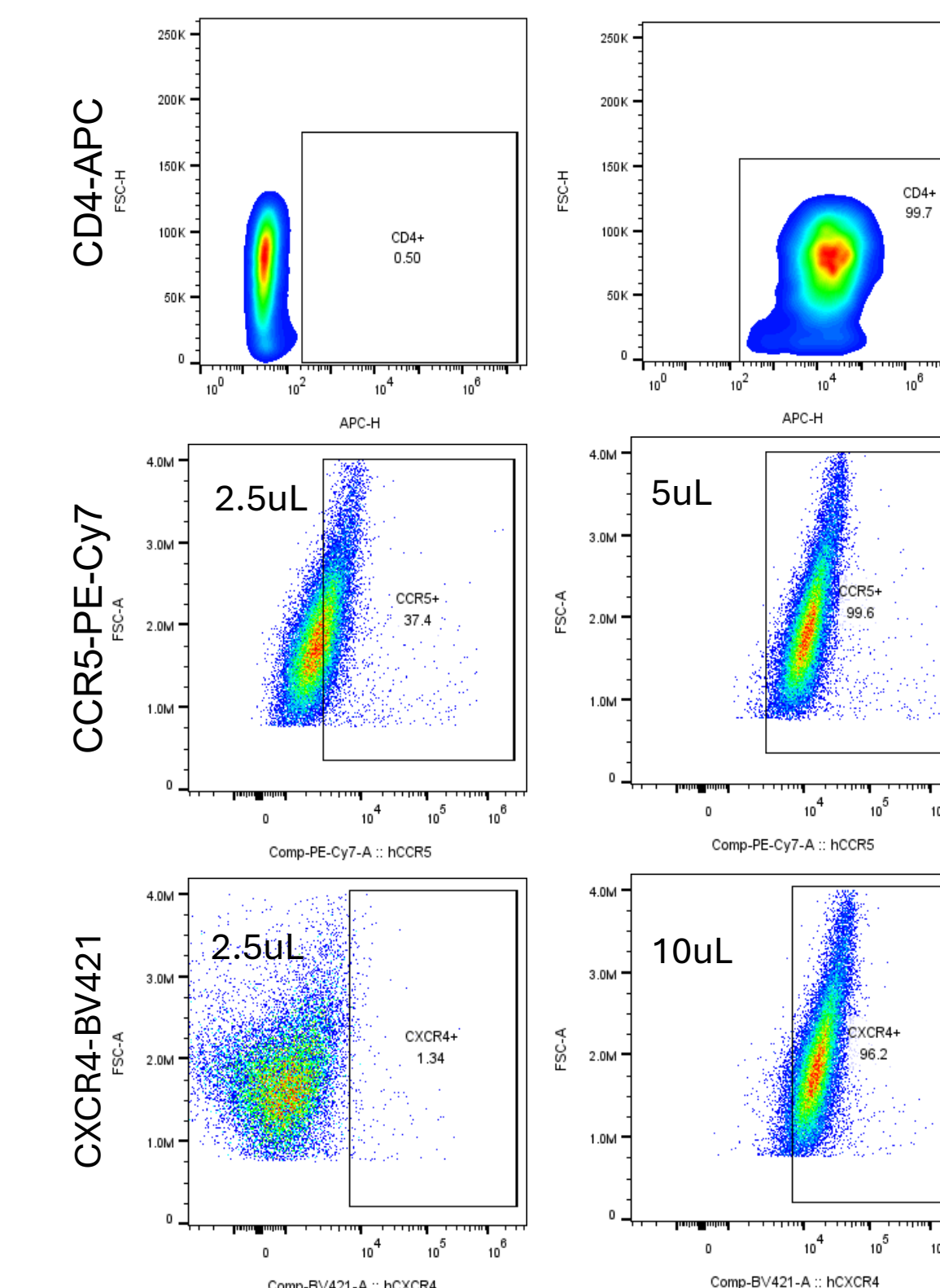


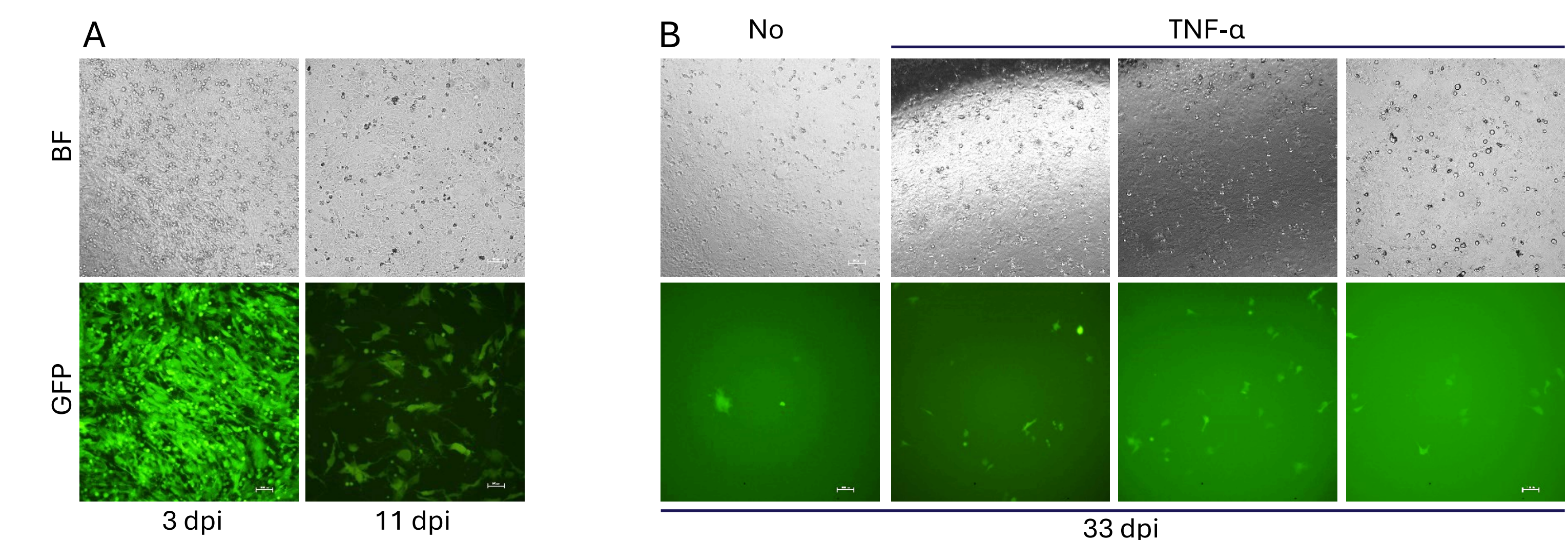
Figure 3. Expression of HIV-1 receptor and co-receptors in HMC3-CD4 clone 9



Proof of Concept: Establishment of HIV Latency in HCM3-CD4 Cells

As a proof of concept to demonstrate that HMC3-CD4 Clone 9 cells can support HIV latency, the cells were infected with a dual-envelope ADA-GFP reporter virus to enhance viral entry. By 3 dpi, ~90% of cells were infected, as determined by GFP expression. However, by 11 dpi, a reduction in GFP expression was observed (Fig. 6A). The cells were passaged for 33 days, after which only 1–2% retained detectable GFP expression. To confirm whether the remaining GFP-negative cells harbored latent HIV, they were treated with TNF- α for 24 hr, followed by GFP monitoring. As shown in Fig. 6B, TNF- α treatment reactivated latent HIV, leading to GFP expression. These data clearly demonstrate that HMC3-CD4 cells can establish HIV latency. Current efforts are focused on generating a single-cell clone of a latent reporter cell line for further studies.

Figure 6. Establishment and Reactivation of Latent HIV in HMC3-CD4 clone 9-6 Cells



Conclusions

- ❖ Human microglia HMC3 cells were successfully constructed for stably expression of HIV CD4 receptor.
- ❖ We demonstrate for the first time that HMC3 cells can support HIV-1 infection, though with low efficiency (~1-2%).
- ❖ Higher CD4 expression minimally enhanced infectivity, as the highest-expressing HMC3-CD4 subclone (9-6) showed only ~5% increased infection, implying additional factors regulate HIV in these cells.
- ❖ HIV latency was successfully established in HMC3-CD4 cells, and current efforts are focused on creating a clonal latent reporter cell line for mechanistic studies and drug screening.

Acknowledgements

This work is supported by an Internal Research and Development (IRAD) program from ATCC. The following reagent was obtained through BEI Resources, NIAID, NIH: Human Immunodeficiency Virus Type 1 (HIV-1) Infectious Molecular Clone, pNL4.3-ADA-GFP-IRES-Nef, HRP-20367.

References

1. Matthew S Dahabieh, Emilie Battivelli and Eric Verdin. Understanding HIV Latency: The Road to an HIV Cure, Annu Rev Med. 2015; Vol. 66:407–421.
2. Yuyang Tang, Antoine Chaillon et al. Brain microglia serve as a persistent HIV reservoir despite durable antiretroviral therapy. Citation Information: J Clin Invest. 2023;133(12):e167417. <https://doi.org/10.1172/JCI167417>
3. Cinzia Russo, Natalia Cappoli et al. The human microglial HMC3 cell line: Where do we stand? A systemic literature review; Journal of Neuroinflammation; 2018 Sep Vol. 10;15:259.