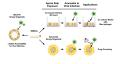
# Apical-Out Airway Organoids and Airway-Immune Co-Cultures Are Suitable Models for Antiviral Compound Screenings and to Study Infectious Diseases

There has been a recent shift away from both in vivo and transformed cell line models and towards advanced primary-derived human in vitro culture systems. Advanced cell models, such as arway organolist and aptihelial-immune cell co-cultures, could serve as powerful platforms to study viela relation and pathogenesis. Despite the advancements in human primary-derivate in vitro models, limitations still remains, including the challenge of accessing the apical surface of epithelia in traditional groanoids with basal side-out polarity, as well as the lack of heterogeneity and paracrine

1. Co-culture model with macrophages.

We developed a workflow to generate comp co-cultures of human bronchial epithelial or (HBECs) and blood-derived macrophages. HBECs

 Apical-out einway organoids for drug screening.
The establishment of an air-liquid interface culture matrix (ECM)-embedded airway organoids, the apical side faces inward. To address these limitations, we

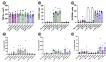


## FIGURE 1. Two Human Airway Model Systems for Viral Infection.

Diagram illustrating two apical-out models for viral infection described in this work: A 2D monolayer model on cell culture inserts for co-culture systems (top), and airway organoids for drug screening (bottom).

Macrophage Culture: Human blood-derived monocytas were isolated using EasySep<sup>TM</sup> Monocyte Isolation Kit and cultured in ImmunoCult<sup>IM</sup>-SF Macrophage Medium (MM), before differentiation to naive (MD-like) macrophage using MD medium (MM - 50 ng/mL macrophage colory-stimulating tactor) for 3 - 4 days. MD site macrophage were turther differentiated for 2 - 1 days into M1 - or M2-like macrophage using M1 medium (MD medium + + 50 ng/mL TMP + 10 ng/mL Experimentation or M2 medium (MD medium + 10 ng/mL L - 4), respectively.

Co-culture and RSV infection: Differentiated air-liquid interface cultures were generated from primary-issue darived HBECs, which were first expanded in PneumaCu<sup>HM</sup>-ExPlus and subsequently differentiated in PneumaCu<sup>HM</sup>-ALI (ALI) medium. After 4 week of differentiation at the air-liquid interface, HBECs in Transvelle inserts were placed on top of differentiated mecophages at the bottom of the lower compartment in either HildFlack, HECL3 in Trainwall® riferra were packed on op a sinstreaman macropragia is in bootin a service control of the service of the se of HRECs for harrier function



### FIGURE 2. HBECs and Macrophages in Co-culture Were Not Significantly Affected by the Cell Culture Medium or Presence of the Other Cell Type.

(A) TER measurements for HBECs were not significantly affected by buture in different microphage media, nor by co-culture with microphages. (B) Flow cytometry analysis revealed that M microphages expressed M1-like marker CD80, and that (C) M2 macrophages expressed the M2 marker CD209 with no significant changes in these markers in ALI the M2 marker CL2car with no argument changes in these measures in the medium or after co-culture. Macrophage phenotype was confirmed by EUSA through release of (D) IL-12 and (E) TNFe for M1-like macrophages, or (F) IL-10 for M2-like macrophages. No significant changes were observed after co-culture. Data represent mean a SEM (n = 4 - 5 donors)

### FIGURE 3. Co-culturing HBECs with M1 Macrophages Produced a Decrease in RSV Infection in HBECs.

(A) Representative images of infected co-cultures. The resulting fluorescence in ALI medium compared to MD or M2 media, indicating that the choice of 5 donors). Scale bar = 1 mm. (B) Quantification of NS1 by oPCR revealed co-culture with M1 macrophages in IMM or M1 medium significantly reduced RSV infection compared to infection in IMM without co-culture. Increased RSV infaction was observed in co-cultures of MD in IMM and co-cultures of M2 in M2 medium, compared to IMM alone, ALI medium reduced the ability of M1 macrophages to reduce infaction. \* = compared to IMM, # = compared to ALJ Imean a SEM, n = 3 - 5 donors). For statistical analysis, ANOVA was performed with Tukey's or Dunnett's post hoc tests as appropriate. \* or # = p < 0.05. \*\* or ## = p < 0.01. \*\*\* or ### = p < 0.001. \*\*\*\* or #### = p < 0.0001

Generation of AOAOs: HBECs were cultured in PneumaCult<sup>IIII</sup>-Ex Plus Medium in 2D, dissociated into single cells, resuspended in PneumaCult<sup>IIII</sup> AOAO Medium, and seeded at 100 cells/microwell of an AggreMell<sup>IIII</sup>400 24-well plate pre-treated with Anti-Adherence Finsing Solution. Seeded cells generated Worder, and swedels as foo committeewale of an egginitive "wood 24-wall place pre-instand with "Unit-infinitements interant good and aggregates after 1 - 6 days in the Aggregate (Well "Place. Aggregates were resultened in the same culture medium and transferred to a new 24-well non-adherent. Taking culture clate for adversarie subprintion culture. Mart 9 - 14 days of subprintion culture. Enrichmal citia were observed at the outer side of the organistic.

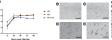
Wrai Infection: Day 15 ADADs were infected with influenza A, influenza B, and rhinovirus A16 at 0.2 MDI and enterovirus D88 at 0.1 MDI. To quantify viral RNA. load, 100 uL of spent medium was sampled at 24-hour intervals for 96 hours and assessed by oPCR. Cultures were topped up with fresh medium after

Antivinal Assay: To datermine the effect of antivinal compounds on the vinal load of enterovinus DB8-infected ACARs, 2.5 µM rupintivir or 5 µM traconazole was added to the cultures at 0 hours post-infection. 100 µL of medium was collected to datect viral RNA load at 24, 48, and 72 hours post-infection.



### FIGURE 4. Apical-out Organoids have Outward Facing Actively Beating Cilia and Express Expected Airway Ciliated Cell Markers Similar to Air-Liquid Interface Cultures,

(A) Representative brightlind image of a day 15 offlorentiated ADACs with cilis on the outer surface. (B) Fundacent immunothermistry analysis of by 15 offlorentiated ADACs with cilis on the outer surface. (B) Fundacent immunothermistry analysis of built of the advance of mining the approximation markers in ADACs with cilis on the outer surface confirming the approximation markers in ADACs with cilis of the approximation and the advance of the advanc Expansion Medium (PreumaCut<sup>IN</sup>-Ex Plus), and at a similar level as 2D differentiated cultures on inserts in ALI medium. Low expression of secretory ineages (i.e. MUCSAC) was detected in AOADs compared to ALI medium. Scale bars = 50 µm.





(A) Successful viral infection and replication were observed in AOADs using epresented as copiessample, log10 of the mean a SD (n = 3) Cytopathogenic effect (CPE) following infection with (B) IAV, (C) IBV, and (D) RV-A16, compared to (E) mode-infacted AOADs at 48 hours obst-infacted. AOADs infacted with IAV or IBV displayed high levels of CPE.



### FIGURE 6. AOAOs are a Suitable Model for Antiviral Drug Screenings.

(A) Quantification of the viral RNA load in enterovirus (EV-DB8)-infected ACA/Da with no addition of drug (black bars), with 5 µM inaconazole (ITZ, orange bars) or 2.5 µM repirit/vir (RUP, blue bars). ITZ treatment significantly inhibited in vitro EV-DB8 replecation primitivy during the tract to tours, writewas the compensation abolished replication (mean a SD, n = 3 donors), "= p < 0.01 and "" = p < 0.001 (#) CPE after 72 hours following inflection of ADADs with EV-D68. Infected ADADs treated with animital drugs showed partial to complete CPE rescue when treated with (C) ITZ and (D) RUP, respectively, Scale bars = 300 um

- · Our work showcases two different models to study viral infection with different characteristics.
- \* A co-culture system that can accommodate assays requiring increased complexity with its capacity to model epithelial and
- . A cost-effective, completely defined, and scalable workflow for AOAD generation that overcomes the limitations of current primary
- Country from HBECs with M1 macrombanes considerably decreased the initial RSV infection, while country with M0 or M2
- \* These results demonstrate that different macrophage subsets play different roles during RSV infection, and these interactions
- ADAOs generated with this ECM-free, defined, and scalable workflow;
- Are susceptible to infection with influenza A. influenza B. rhinovirus A16, and enterovirus D68.
- . Successfully modeled the antiviral effect of rupintrivir and itraconazole in viral RNA load and CPE following enterovirus D68.



