

Forskolin-Induced Swelling of Human Intestinal Organoids Grown in IntestiCult™

Cystic fibrosis (CF) is a genetic disease caused by defects in the gene encoding the cystic fibrosis transmembrane conductance regulator (CFTR), a chloride anion channel expressed in multiple epithelial tissues, including the intestinal epithelium. Defects in CFTR cause aberrant ion transport and mucus build-up in various organs, leading to difficulty breathing, recurring lung infections, and digestive issues. The recent characterization of intestinal organoid cultures has provided a new tool for studying CFTR function in vitro.¹ Intestinal organoids can be efficiently established from colorectal biopsies, expanded, and maintained in long-term in vitro culture. These organoids maintain the genotype and phenotype of the parental tissue and retain the expected CFTR function when cultured in vitro.

Exposing intestinal organoids to forskolin causes the cells to rapidly increase their levels of cyclic adenosine monophosphate (cAMP), resulting in the opening of the CFTR channel.^{1,2} As chloride ions move through the channel and into the lumen, the organoids increase in size due to water following by osmosis, allowing for the assay of CFTR activity through organoid swelling. Because the swelling is CFTR-dependent, wild-type organoids increase in size while those carrying a CFTR mutation, or treated with a CFTR inhibitor, do not.

Below we describe a protocol for forskolin-induced swelling of human intestinal organoids derived from healthy individuals and cultured in IntestiCult™ Organoid Growth Medium (Human) (#06010). For complete culturing instructions, use this document in conjunction with the relevant Product Information Sheet (Document #DX21423).

Materials

Product Name	Catalog #
Forskolin	72112
CFTR _{INH} 172	e.g. Cayman 15545
GlyH-101	e.g. Cayman 15772
DMSO	-
Costar® 24 Well Flat-Bottom Plate, Tissue Culture-Treated	38017
Krebs-Ringer Solution, Bicarbonate-Buffered	e.g. Alfa Aesar J67591

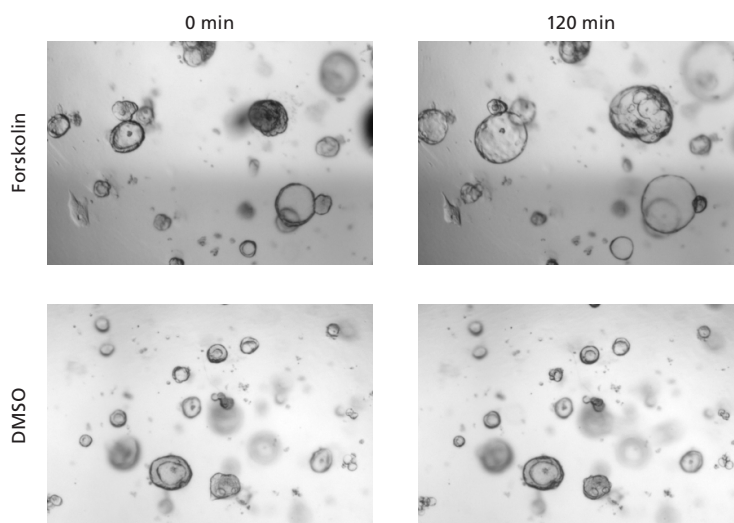


Figure 1. Forskolin-Induced Swelling of Intestinal Organoids Cultured in IntestiCult™ Organoid Growth Medium (Human)

The lumen of intestinal organoids became enlarged in response to treatment with 10 μ M of forskolin after 120 minutes. This increase in size was not observed in intestinal organoids treated with a DMSO control.

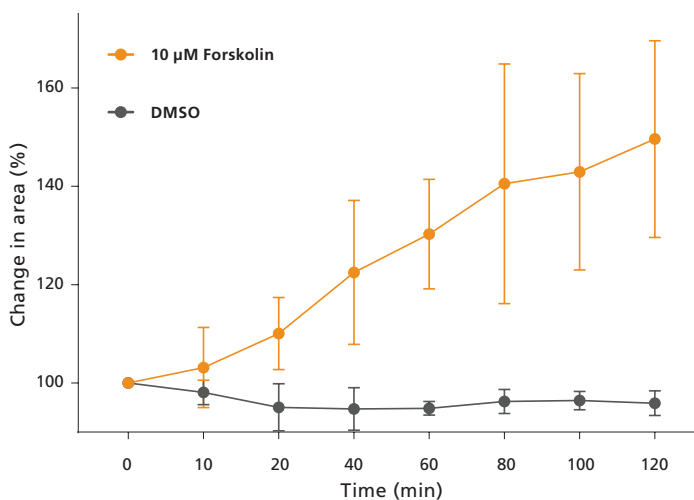


Figure 2. Quantification of Forskolin-Induced Swelling of Intestinal Organoids

The swelling of intestinal organoids was quantified by brightfield microscopy. The percentage change in area at later time points was calculated relative to the average area of the organoids observed at time = 0. At each time point, the whole well was imaged and the average from three technical replicates was calculated. Data is representative of one normal colonic donor at passage 3.

Forskolin-Induced Swelling of Intestinal Organoids

1. Establish organoid cultures as described in the Product Information Sheet (Document #DX21423). Grow organoids for 3 - 5 days after passage in IntestiCult™ Organoid Growth Medium (Human) (Catalog #06010) prior to assessing forskolin-induced swelling.
2. Remove media from each well and replace with an equal volume of Krebs-Ringer Bicarbonate (KBR) buffer. Allow the buffer to equilibrate with the Corning® Matrigel® domes for 30 minutes at 37°C.

NOTE: Forskolin-induced swelling can be performed and assayed in the culture medium, however, swelling is more pronounced if using KBR buffer.
3. OPTIONAL: If using CFTR inhibitors as an additional control (e.g. CFTR_{INH}172 or GlyH-101), add the appropriate amount of inhibitor to each well as necessary (50 µM for CFTR_{INH}172 or GlyH-101) and incubate for 3 hours at 37°C. If inhibitors are not being used, proceed directly to step 4.
4. Capture time = 0 images before proceeding as organoid swelling may begin immediately after adding forskolin.
5. Remove the buffer from all wells being assayed and replace with fresh buffer, inhibitors (if applicable), and forskolin to a final concentration of 10 µM forskolin.
6. Image each well at regular intervals until the swelling in uninhibited wells ceases (usually 80 - 120 minutes).
7. To quantify forskolin-induced swelling, brightfield images at each time point are analyzed using ImageJ for average organoid area compared to the time = 0 images.

Related Products

Description	Quantity	Catalog #
IntestiCult™ Organoid Growth Medium (Human)	100 mL	06010
IntestiCult™ Organoid Growth Medium (Mouse)	100 mL	06005

References

1. Dekkers, JF et. al (2012). A functional CFTR assay using primary cystic fibrosis intestinal organoids. Nat. Med. 19(7): 939–47.
2. Boj, SF et. al (2017). Forskolin-induced swelling in intestinal organoids: An in vitro assay for assessing drug response in cystic fibrosis patients. Journal of Visualized Experiments. 120(e55159).

Copyright © 2020 by STEMCELL Technologies Inc. All rights reserved including graphics and images. STEMCELL Technologies & Design, STEMCELL Shield Design, Scientists Helping Scientists, and IntestiCult are trademarks of STEMCELL Technologies Canada Inc. Corning, Costar, and Matrigel are registered trademarks of Corning Incorporated. All other trademarks are the property of their respective holders. This product was developed under a license to intellectual property owned by The Hubrecht Organoid Technology Foundation ("the HUB"). This product is sold for research use only. Purchase of this product does not include the right to use this product for growing organoids for drug screening aiming for commercial gain or for other commercial purposes. Purchasers wishing to use the product for purposes other than research use should contact the HUB to obtain a further license (an "Organoid-Growth License"). Purchasers may apply for an Organoid-Growth License from the HUB and such license will not be unreasonably withheld by the HUB. While STEMCELL has made all reasonable efforts to ensure that the information provided by STEMCELL and its suppliers is correct, it makes no warranties or representations as to the accuracy or completeness of such information.

STEMCELL TECHNOLOGIES INC.'S QUALITY MANAGEMENT SYSTEM IS CERTIFIED TO ISO 13485. PRODUCTS ARE FOR RESEARCH USE ONLY AND NOT INTENDED FOR HUMAN OR ANIMAL DIAGNOSTIC OR THERAPEUTIC USES UNLESS OTHERWISE STATED.