

Optimization of tumor tissue enzymatic dissociation and initial culture conditions for rapid and viable tumoroid generation

Logan Wilson, Colin Paul, Anthony Chatman, Chris Yankaskas, Pradip Shahi Thakuri, Matt Dallas, and David Kuninger
Thermo Fisher Scientific, Frederick, MD, USA 21704

Key Takeaways

- OncoPro™ Tumoroid Culture Medium is a standardized, serum-free and conditioned medium-free media system for patient-derived tumoroid culture
- Optimization of tumor dissociation conditions balances viable tumor cell yield with tumoroid formation in culture
- Addition of growth factors during initial derivation may improve long-term tumoroid establishment efficiency

Introduction

Several studies have indicated that patient-derived cancer models (tumoroids or cancer organoids) recapitulate key genotypic and phenotypic features of patient tumors, making them potentially valuable tools for basic and translational cancer research. OncoPro™ Tumoroid Culture Medium is a serum-free, conditioned medium-free cell culture medium designed to culture tumoroids derived from multiple cancer indications. Here, we highlight studies aiming to improve tumoroid formation during initial culture establishment, focusing on the development of enzymatic dissociation cocktails that maximize tumoroid formation in culture and results of screening experiments designed to improve tumoroid formation during initial tumoroid passaging.

Materials and Methods

Surgical resection samples were placed in supplemented Hibernate™-A Medium containing antibiotics and shipped overnight at 4° C to a central processing lab. Tissues were dissociated via enzymatic and mechanical dissociation using various enzyme cocktails, counted, and plated in OncoPro™ Tumoroid Culture Medium (in the presence of indication-specific factors when applicable). Cultures were analyzed for metabolic activity (using PrestoBlue™ HS Cell Viability Reagent) and tumoroid area (by imaging on an EVOS™ M7000 Imaging System) after 7 days in culture. The workflow is illustrated schematically in Figure 1. In separate experiments, dissociated tumor cells from resections or commercial sources were cultured in OncoPro™ Tumoroid Culture Medium supplemented with various factors, and tumoroid size and cell count were tracked over initial passage(s).

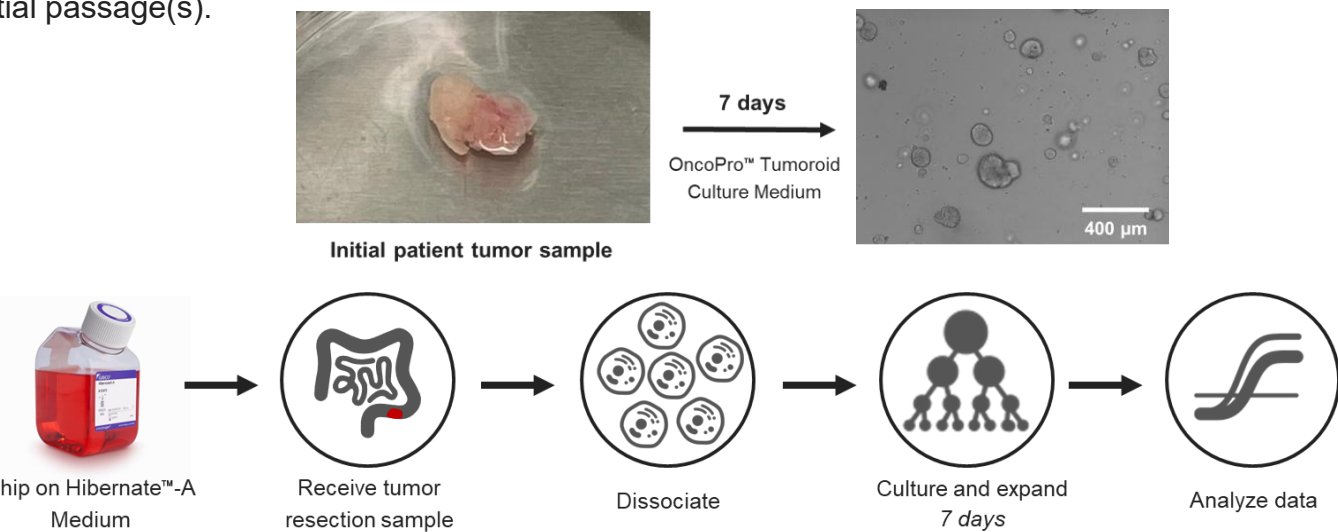


Figure 1. Experimental workflow for generation of tumoroid from cancer tissue samples. Surgical resection samples were shipped overnight in supplemented Hibernate-A medium, dissociated, and plated in OncoPro Tumoroid Culture Medium. After 7 days of culture, the resulting tumoroids were analyzed.

Results

Transport medium design for retained cell viability during tissue shipment and storage

Tissue transport buffer was designed to promote cell survival in CO₂-independent conditions. Selected samples were divided into pieces upon receipt and stored in the tissue transport buffer at 4°C for various intervals prior to enzymatic dissociation and quantification of cell number and viability. Samples retained similar viability during storage in transport buffer (Figure 2).

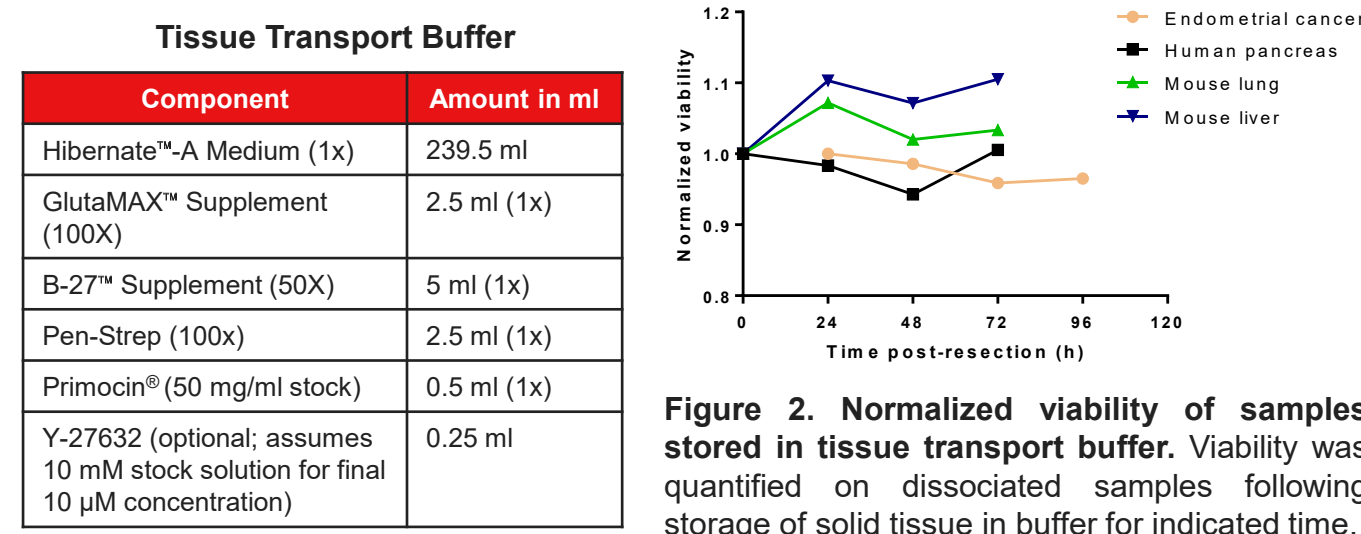


Figure 2. Normalized viability of samples stored in tissue transport buffer. Viability was quantified on dissociated samples following storage of solid tissue in buffer for indicated time.

Tissue dissociation workflow

Following transport, tissues were placed in tissue processing buffer containing enzymes that promote viable cell release. Specifically, received tissues were minced, mixed well, and divided for dissociation in multiple enzyme mixtures as dictated by a design-of-experiment approach to study a range of components and concentrations in a donor-blocked format (Figure 3). Dissociation data pooled by donor across conditions indicated effective dissociation, with a wide range in post-dissociation yield based on cancer indication. When an identical number of viable cells were seeded following dissociation in distinct enzyme mixtures, differential tumoroid formation was observed. Further analysis of enzymatic dissociation cocktails leading to high yields and tumoroid formation in culture are discussed in the next section.

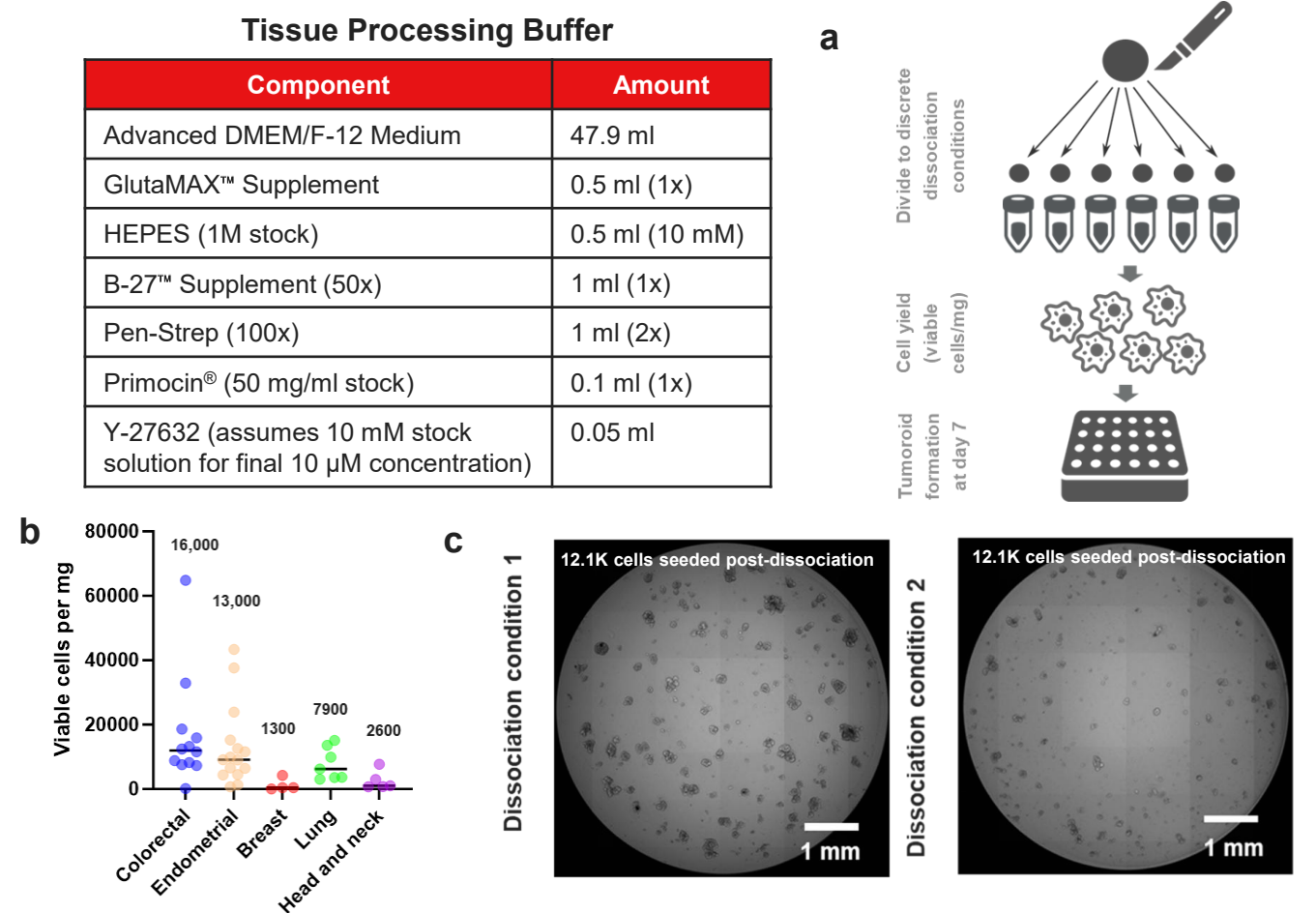


Figure 3. Cell yields and viabilities following dissociation of solid tumor resection samples. (a) Tissue dissociation workflow. (b) Viable cells per mg tissue after tissue dissociation (pooled across conditions for a given donor). Mean viable cells per mg is displayed for each indication. Each point represents one donor. (c) Representative images of tumoroid formation following dissociation of donor-matched tissue with differing enzymes.

Optimization of tissue dissociation cocktails

Tumor tissue dissociation cocktails were determined based on analysis of viable cell yield and health in culture for multiple donors per cancer indication. Example data for colorectal cancer samples are shown below (Figure 4). Final enzymatic cocktail formulations for four cancer indications are shown.

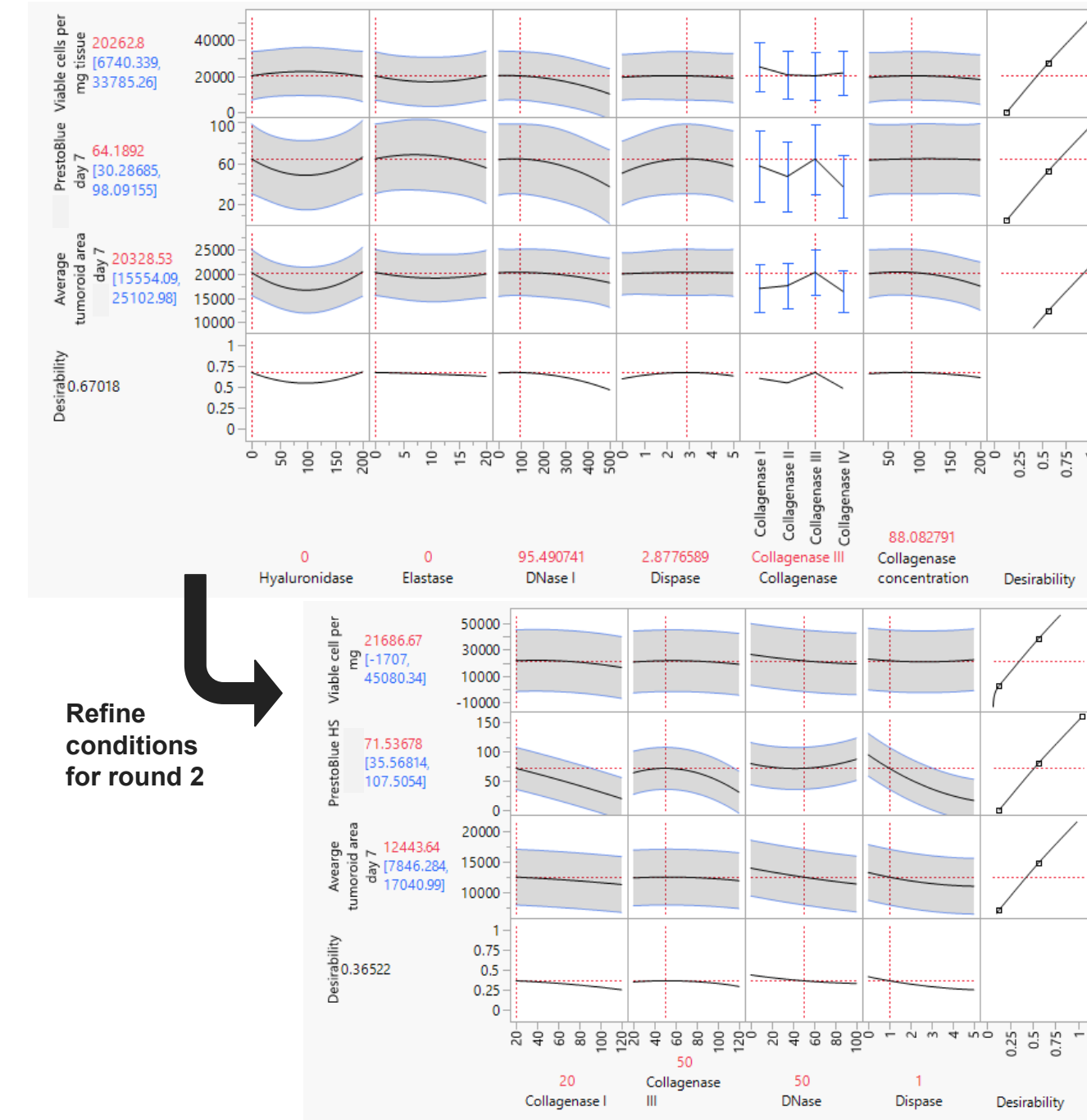


Figure 4. Effect of enzyme concentration on tissue dissociation and initial colorectal tumoroid culture. Impact of enzymes on viable cell yield post-dissociation and tumoroid formation following seven days of culture. Five donors were used for each optimization round.

Colorectal cancer		Lung cancer	
Component	Amount	Component	Amount
Tissue Processing Buffer	2.5 ml	Tissue Processing Buffer	2.5 ml
Collagenase Type I	20 U/ml	Collagenase Type I	80 U/ml
Collagenase Type III	50 U/ml	Hyaluronidase	50 U/ml
Dispase	1 U/ml	Dispase	1 U/ml
DNase I	50 U/ml	DNase I	100 U/ml

Breast cancer		Endometrial cancer	
Component	Amount	Component	Amount
Tissue Processing Buffer	2.5 ml	Tissue Processing Buffer	2.5 ml
Collagenase Type III	150 U/ml	Collagenase Type III	5 U/ml
Dispase	1 U/ml	Collagenase Type IV	5 U/ml
DNase I	50 U/ml	Dispase	7 U/ml
		DNase I	50 U/ml

Screening of factors to promote tumoroid expansion during initial culture

Viable cryopreserved dissociated tumor cell (DTC) samples and dissociated breast cancer samples were cultured in OncoPro medium supplemented with additional factors in a designed experiment. Cells were imaged over time for analysis of tumoroid area, and tumoroids in each condition were dissociated and counted. The influence of each input supplement was modeled to identify factors that potentially increase tumoroid formation and cell number (Figure 5).

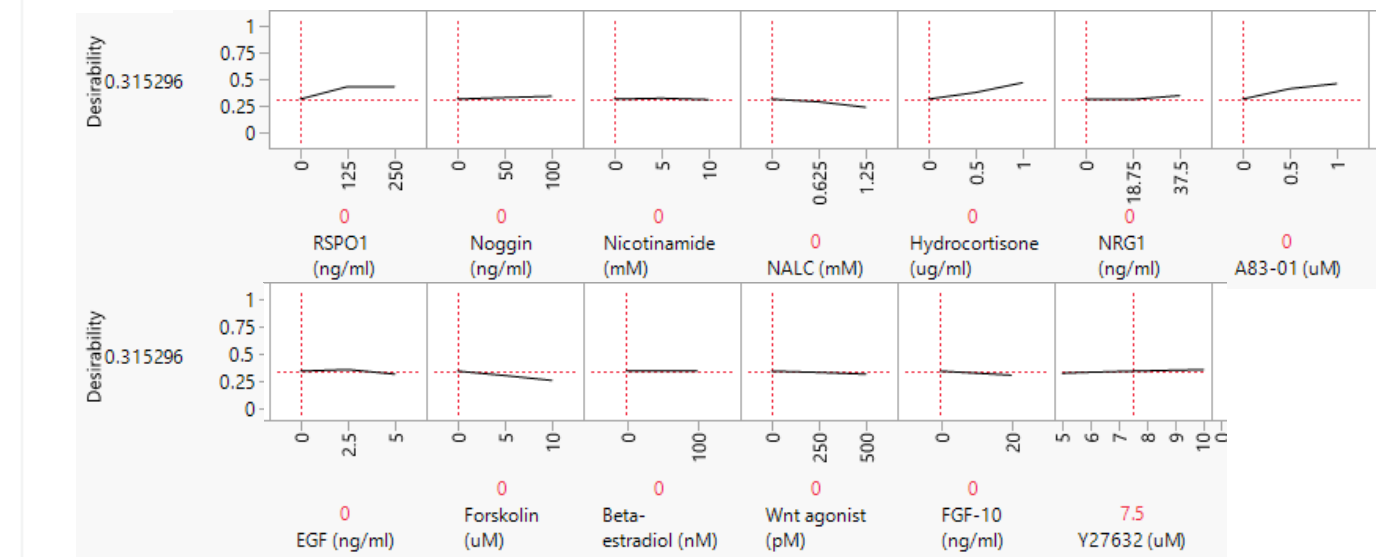


Figure 5. Effects of supplements to OncoPro™ Tumoroid Culture medium. Desirability plots were generated from data on tumoroid area and cell count during initial breast cancer cell culture. Results are consolidated across six donors.

Following screening for factors promoting viable tumoroid cells during initial culture, a prototype formulation was generated and tested on a fresh breast cancer tissue sample, on which cell counts were tracked over multiple passages (Figure 6).

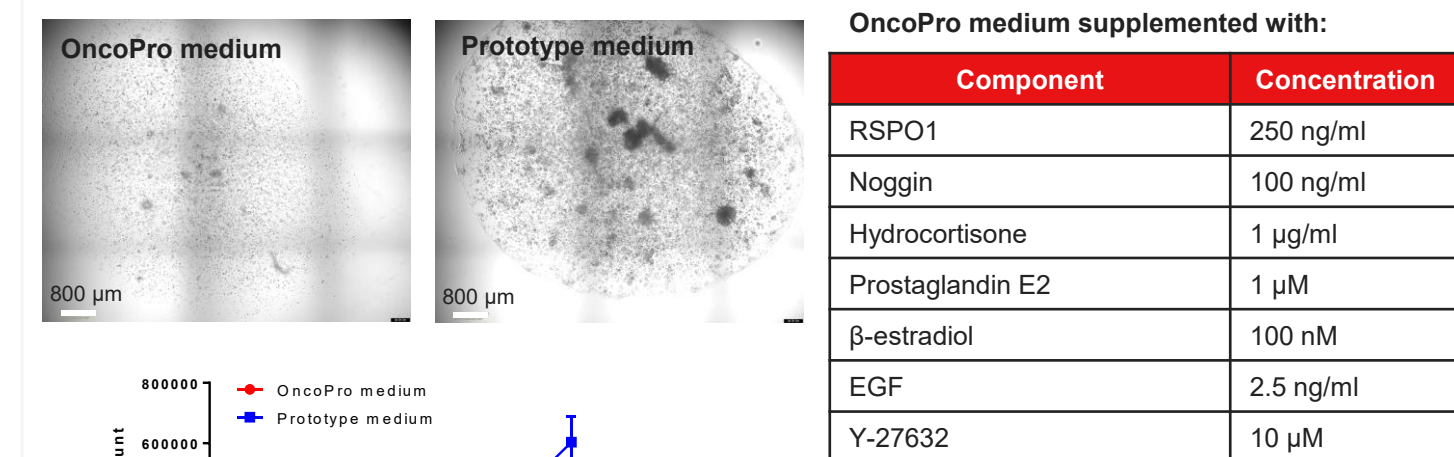


Figure 6. Breast tumoroid expansion in prototype derivation medium. Representative images after 13 days of culture and cell counts over time following culture in OncoPro medium for breast cancer (includes 10 ng/ml heat stable FGF-10 and 10 nM β-estradiol) or supplemented OncoPro medium.

Trademarks/licensing

© 2026 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others. **For Research Use Only.** Not for use in diagnostic procedures.

Science at a scan
Scan the QR code on the right with your mobile device to download this and many more scientific posters.