

Product datasheet

Alginate Hydrogel Kit for 3D Cell Culture ab241011

[1 Image](#)

Overview

Product name	Alginate Hydrogel Kit for 3D Cell Culture
Assay type	Cell-based (qualitative)
Product overview	Alginate Hydrogel Kit for 3D Cell Culture (ab241011) offers 3D cell culture matrices to meet the needs and requirements of various research fields. This kits provide a standardized, yet user friendly and adaptable to high-throughput strategy for setting up spheroid formations, 3D cell cultures and pharmacological studies.

Properties

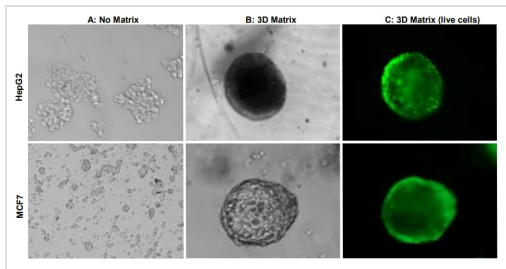
Storage instructions Store at -20°C. Please refer to protocols.

Components	100 tests
Alginate Hydrogel Matrix	1 x 5ml
Cross-linking Solution	1 x 25ml
Wash Buffer	1 x 100ml

Relevance

Three dimensional (3D) cell cultures are artificially-created environments in which cells are permitted to grow or interact with their surroundings in a 3D fashion. 3D cell cultures improve the function, differentiation and viability of cells and recapitulate in vivo microenvironment compared to conventional 2D cell cultures. 3D matrices provide a physiologically relevant screening platform, by mimicking the in vivo responses, for many cell types including cancer and stem cells in developmental morphogenesis, pharmacology, drug metabolism and drug toxicity studies.

Images



HepG2 and MCF7 cells in No Matrix (A) and 3D Alginate Hydrogel Matrix (B).

HepG2 and MCF7 cells in No Matrix (A) and 3D Alginate Hydrogel Matrix (B). Cells were cultured in Alginate Hydrogel Matrix for 21 days, and successfully formed spheroids. Media was changed every 2-3 days as per protocol. The Calcein AM staining (C) indicates that cell viability is not affected while culturing in matrix for a long period of time. Note: Calcein AM is not included in the kit.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors