

Product datasheet

3D Cell Culture Clearing Kit ab243299

2 Images

Overview

Product name 3D Cell Culture Clearing Kit

Product overview 3D Cell Culture Clearing Kit ab243299 enables the easy set-up of rapid 3D cell culture clearing for 3D imaging using simple protocols and standard laboratory equipment.

This kit can be used with immunostaining, fluorescent proteins and chemical dyes. Clearing is reversible so that you can treat your sample for conventional H&E or IHC staining after 3D imaging.

To help you get started with clearing 3D cell cultures, we have [clearing validated antibodies](#). Validated antibodies are for key markers, such as NeuN, GFAP, Iba1, etc, and are tested for [tissue clearing](#) with 1 mm mouse brain sections, and for 3D cell culture clearing with neuronal spheroids and hepatic spheroids (depending on the marker).

If you want to set up clearing of 3D cell cultures on a sample type that we haven't tested or to use your own antibodies, please consult the protocol book for this kit.

To learn more about tissue clearing and clearing of 3D cell cultures, please review our [guide to tissue clearing and clearing 3D cell cultures](#).

[Clearing reagents and companion products](#) are also available as individual products.

These clearing reagents were developed with Visikol Inc, who developed the Visikol® HISTO™ clearing technology which is the basis of our reagents.

Tested applications **Suitable for:** Histology

Properties

Storage instructions Please refer to protocols.

Components	30 ml	15 ml
Tissue Clearing Antibody Buffer	1 x 30ml	1 x 15ml
Tissue Clearing Blocking Buffer	1 x 30ml	1 x 15ml

Components	30 ml	15 ml
Tissue Clearing Penetration Buffer	1 x 30ml	1 x 15ml
3D Cell Culture Clearing Reagent	1 x 30ml	1 x 15ml
Tissue Clearing Washing Buffer 10X	1 x 70ml	1 x 35ml

Relevance

The goal for tissue clearing is making large, fixed biological samples transparent. “Large” in this case means thick sections of tissue, whole organoids, entire organs or even entire young rats, ranging in thickness from around 100 µm to several centimeters. Typically, samples of this size are not transparent, and are therefore difficult to analyze using visible wavelengths of light with a microscope. Tissue clearing involves a series of chemical steps that render a large sample transparent.

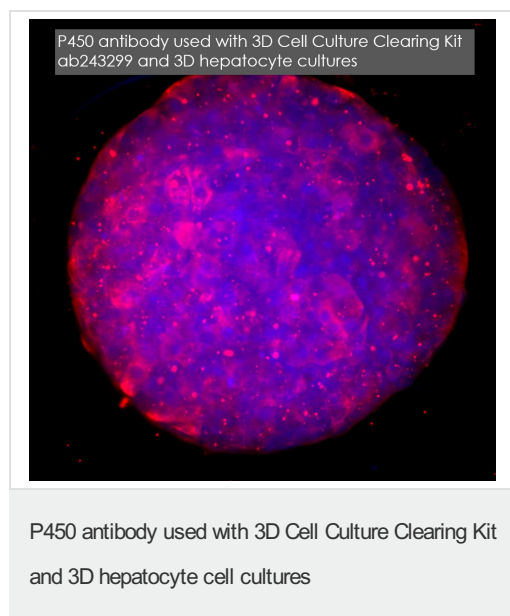
Applications

Our [Abpromise guarantee](#) covers the use of **ab243299** in the following tested applications.

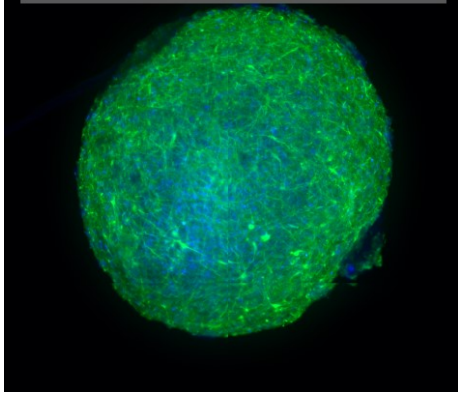
The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Histology		Use at an assay dependent concentration.

Images



GFAP antibody with 3D Cell Culture Clearing Kit
ab243299 and neuronal spheroid cell culture



GFAP antibody

ab7260

with 3D Cell Culture Clearing Kit and 3D neuronal
spheroid cell cultures

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