

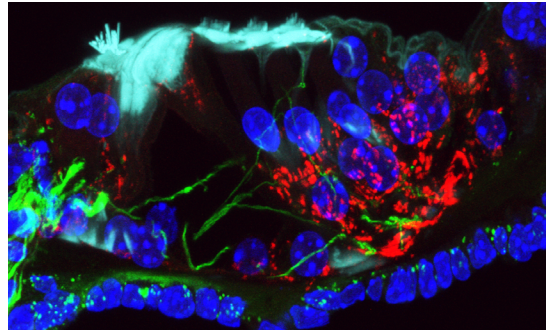
VECTASHIELD® Antifade Mounting Medium with DAPI

Cat. No.: H-1200

Concentration of DAPI: 1.5 µg/ml

Storage: Refrigerate

Refractive Index: 1.45



Adult organ of Corti labelled with anti-beta3-tubulin (green), phalloidin (cyan), anti-connexin30 (red) and DAPI (blue). Image provided by Dan Jagger, Ear Institute, University College London. Mounted in VECTASHIELD Antifade Mounting Medium with DAPI.

DESCRIPTION

VECTASHIELD Antifade Mounting Medium is a unique, stable formula for preserving fluorescence. VECTASHIELD Antifade Mounting Medium prevents rapid photobleaching of fluorescent proteins and fluorescent dyes (see website for compatibility).

Each lot is tested for photobleaching retardance with qualitative and quantitative methods.

This formulation of VECTASHIELD Antifade Mounting Medium contains 4',6-diamidino-2-phenylindole (DAPI), which fluoresces when bound to DNA. DAPI excites at about 360 nm and emits at about 460 nm when bound to DNA, producing a blue fluorescence, and is ideal as a nuclear or chromosomal counterstain.

APPLICATION

To mount tissues or cells on a slide dispense VECTASHIELD Mounting Medium onto the specimen using a disposable pipet or pipet tip. Small drop volumes of approximately 25 µl (per 22 mm x 22 mm coverslip) are recommended. Coverslip the specimen and allow VECTASHIELD Mounting Medium to disperse over the entire section. Tightly seal the bottle with the screw cap after use and store in an upright position.

VECTASHIELD Antifade Mounting Medium does not solidify, but remains a liquid on the slide and can be stored without sealing. If desired, coverslips can be permanently sealed around the perimeter with nail polish or a plastic sealant. Mounted slides should be stored at 4 °C protected from light.

Note: DAPI is a mutagen and appropriate precautions should be taken when using it.

VECTASHIELD Antifade Mounting Media Products:

Product	Counterstain	Cat. No.	Unit Size	Hardening
VECTASHIELD Antifade Mounting Medium	none	H-1000	10 ml	no
	DAPI	H-1200	10 ml	no
	PI	H-1300	10 ml	no
VECTASHIELD HardSet™ Antifade Mounting Medium	none	H-1400	10 ml	yes
	DAPI	H-1500	10 ml	yes
	TRITC-Phalloidin	H-1600	10 ml	yes
VECTASHIELD Vibrance™ Antifade Mounting Medium	none	H-1700	2 ml, 10 ml	yes
	DAPI	H-1800	2 ml, 10 ml	yes