

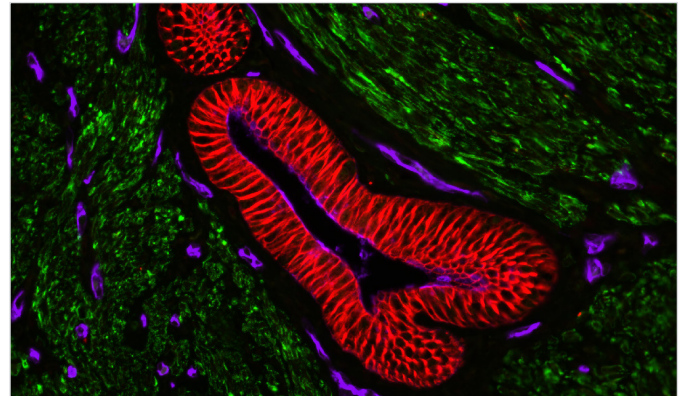
VECTASHIELD Vibrance®

Antifade Mounting Medium

Cat. No.	H-1700
Storage	Refrigerate
Refractive Index	1.38 (initial) 1.47 (cured)
Description	VECTASHIELD Vibrance Antifade Mounting Medium prevents rapid photobleaching of fluorescent proteins and fluorescent dyes on cell and tissue section preparations. VECTASHIELD Vibrance Antifade Mounting Medium is supplied as an aqueous, ready-to-use formulation which will cure (set) over several hours at room temperature to immobilize the coverslip. No further sealing of the coverslip is required.

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

VECTASHIELD Vibrance Antifade Mounting Medium is compatible with the majority of commercially available fluorescent dyes. The antifade properties of the medium are effective across the spectrum, including far red channels.



Human Uterine section (FFPE): Stained for Desmin (green) and Cytokeratin (red) using VectaFluor™ Duet Double Labeling Kit (DK-8828), and vasculature using DyLight™ 649 UEA I lectin (purple). Mounted in VECTASHIELD Vibrance Antifade Mounting Medium.

Application

To mount tissue sections or cells, prepare the slide by removing excess liquid. Keep the specimen moist and do not dry. Use a disposable pipet or pipet tip to dispense VECTASHIELD Vibrance Antifade Mounting Medium onto the specimen and coverslip. Small drop volumes of approximately 25 µl (per 22 mm x 22 mm coverslip) are recommended. For specimens mounted on coverslips, dispense mounting medium onto a clean slide and slowly lower the coverslip onto the medium.

Place mounted samples on a flat, dry surface in the dark and allow the medium to cure. Slides can be viewed 1 hour after mounting, however optimal antifade performance is achieved after 2 hours. The refractive index will reach 1.47 when fully cured (24 hours). For long-term storage, store slides at room temperature or refrigerate.

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® PLUS Antifade Mounting Medium	none	H-1900	2 ml, 10 ml	no
	DAPI	H-2000	2 ml, 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

For research use only. Not intended for animal or human therapeutic or diagnostic use.

©2022 Vector Laboratories. All rights reserved.

HardSet, VectaFLuor, VECTASHIELD, and VECTASHIELD Vibrance are trademarks of Vector Laboratories, Inc.

DyLight is a trademark of Pierce Biotechnology, Inc.

H-1700.LBL-02357.Rev01