

eBioscience™ Phalloidin eFluor™ 520

Catalog Number: 59-6559

RUO: For Research Use Only. Not for use in diagnostic procedures.

Product Information

Contents: eBioscience™ Phalloidin eFluor™ 520

REF

Catalog Number: 59-6559

Concentration: 1000X



LOT



Formulation: Frozen liquid, 300 test size

Temperature Limitation: Store at less than or equal to -20°C. Protect from light and moisture.

Batch Code: Refer to vial

Use By: Refer to vial

Description

Phalloidin is a peptide toxin isolated from the *Amanita phalloides* mushroom. Phalloidin selectively binds F-actin and functions to stabilize actin polymers by inhibiting the dissociation of actin monomers from the filament ends. Because F-actin polymers compose the cytoskeleton of almost all eukaryotic cells, phalloidin conjugates can be useful probes for visualizing the structure of live or fixed cells using microscopy. Phalloidin eFluor® 520 can be used to label cells of multiple species. Its excitation and emission is similar to Alexa Fluor 488 and FITC.

Phalloidin eFluor® 520 is supplied as a frozen liquid. To prepare the 1000X stock solution, thaw the vial, add 30 µL high-quality anhydrous DMSO, mix and briefly spin down. The resulting solution (phalloidin DMSO stock) should be aliquoted, protected from light and moisture, and stored at -20°C. To prepare a 1X working solution of phalloidin conjugate, add 1 µL of 1000X phalloidin DMSO stock solution to 1 mL of PBS containing 1% BSA.

Applications Reported

Phalloidin eFluor® 520 has been reported for use in microscopy, immunohistochemical staining, immunocytochemistry, and cell labeling.

Applications Tested

Phalloidin eFluor® 520 has been tested by staining of fixed and permeabilized cells and can be used at 1X for 20-45 minutes at room temperature. Fixation of cells with methanol-containing fixatives prior to phalloidin staining is not recommended. Optimal staining time may be cell type dependent, it is recommended that the incubation time be determined by each investigator for optimal performance in the assay of interest.

eFluor® 520 is a replacement for Alexa Fluor® 488.

Filter Recommendation: When using this eFluor® 520 phalloidin conjugate, we recommend a filter that will capture the 520 nm emission wavelength (for example, Excitation 475/40, 505LP, Emission 535/45). A standard Alexa Fluor® 488 or FITC filter is acceptable.

Special Notes

Staining with Phalloidin eFluor® 520 may be performed at the same time as fluorophore-conjugated antibodies. For best results visualize phalloidin-stained cells within 1 hour of staining.

References

Berrout L, Isokawa M. Ghrelin promotes reorganization of dendritic spines in cultured rat hippocampal slices. *Neurosci Lett.* 2012 May 16;516(2):280-4.

Saraceno GE, Ayala MV, Badorrey MS, Holubiec M, Romero JI, Galeano P, Barreto G, Giraldez-Alvarez LD, Kölliker-Fres R, Coirini H, Capani F. Effects of perinatal asphyxia on rat striatal cytoskeleton. *Synapse.* 2012 Jan;66(1):9-19.

Volin MV, Huynh N, Klosowska K, Reyes RD, Woods JM. Fractalkine-Induced Endothelial Cell Migration Requires MAP Kinase Signaling. *Pathobiology.* 2010 February; 77(1): 7-16.

Related Products

00-4958 Fluoromount-G™

Not for further distribution without written consent.

Copyright © 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

Tel: 888.999.1371 or 858.642.2058 • Fax: 858.642.2046 • thermofisher.com/ebioscience • info@ebioscience.com

eBioscience™ Phalloidin eFluor™ 520

Catalog Number: 59-6559

RUO: For Research Use Only. Not for use in diagnostic procedures.

00-4959 Fluoromount-G™, with DAPI

65-0880 eBioscience™ DRAQ5™

Not for further distribution without written consent.

Copyright © 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

Tel: 888.999.1371 or 858.642.2058 • Fax: 858.642.2046 • thermofisher.com/ebioscience • info@ebioscience.com