

Technical Bulletin

ANTI-FLAG® M2-Cy3 Antibody, Mouse Monoclonal

Clone M2, purified immunoglobulin, buffered aqueous solution

A9594

Product Description

Monoclonal ANTI-FLAG® M2-Cy3 antibody is a covalent conjugate of a purified IgG1 monoclonal ANTI-FLAG® M2 antibody, which has been isolated from a mouse cell culture,¹ to the cyanine dye Cy3™. The antibody conjugate binds to FLAG® fusion proteins. It will recognize the FLAG® sequence at the N-terminus, Met-N-terminus, or C-terminus of FLAG® fusion proteins. This conjugate is useful for identification of FLAG® fusion proteins by common immunological procedures.

Monoclonal ANTI-FLAG® M2-Cy3™ may be used for the detection of FLAG® fusion protein in applications such as fluorescent immunocytochemistry. Cy3™ conjugates are well suited for dual-labeling experiments.^{2,3} Several theses⁴⁻¹⁰ and dissertations¹¹⁻¹⁶ cite use of this product in their research protocols.

Reagent

This product is supplied as a solution in 10 mM sodium phosphate, pH 7.4, with 150 mM NaCl, 1% bovine serum albumin, and 15 mM sodium azide.

Conjugate protein concentration: ~1 mg/mL (exact value on lot-specific Certificate of Analysis)

F/P molar ratio: 3.0-6.0

Specificity: This antibody has been found to detect FLAG-BAP™ fusion protein in transfected COS cells.

Precautions and Disclaimer

For Research use only. Not for drug, household, or other uses. Because of the sodium azide content, a Safety Data Sheet for this product has been sent to the attention of the safety officer of your institution. Consult the Safety Data Sheet for information regarding hazardous and safe handling practices.

Storage/Stability

Store undiluted antibody at -20 °C in working aliquots. Repeated freezing and thawing is **not** recommended.

Preparation Instructions

Dilute the antibody to the recommended working dilution in Tris-Buffered Saline [TBS; 0.05 M Tris (pH 7.4) with 0.15 M NaCl].

Suggested working dilutions:

Immunocytochemistry: an antibody concentration of 1-10 µg/mL may be used.

Note: To obtain best results, it is recommended that each individual user determine the optimal working dilution by titration assay.

Procedure

Procedure for Direct Immunofluorescent Staining of Mammalian Cells

1. Wash cells grown in a culture dish or on a slide with TBS twice.
2. Fix cells with a freshly prepared mixture of methanol:acetone (1:1) for 1 minute at room temperature.
3. Wash cells with TBS four times.
4. Incubate cells with the antibody at 10 µg/mL in TBS at room temperature for 1 hour.
5. Wash cells with TBS twice.
6. Examine using a fluorescence microscope with appropriate configuration for Cy3™. Cy3™ has an absorption maximum at approximately 550 nm with an emission maximum at 570 nm.

References

1. Brizzard, B.L. *et al.*, *BioTechniques*, **16(4)**, 730-735 (1994).
2. Sargent, P.B., *Neuroimage*, **1(4)**, 288-295 (1994).
3. Southwick, P.L. *et al.*, *Cytometry*, **11(3)**, 418-430 (1990).
4. Chun, Sophia Weng Shum, "Overexpression of the transcription factor ZFP60 in a multipotential mesenchymal cell line and its effects on gene expression and cell differentiation potential". University of Toronto, M.Sc. thesis, p. 110 (2008).
5. O'Connell, Marie, "Development of antibodies to human transient receptor potential vanilloid 1 for future targeting of therapeutics to sensory neurons". Dublin City University, M.Sc. thesis, p. 28 (January 2014).
6. Husić, Mia, "Binding, activation and cellular actions of teneurin C-terminal associated peptide (TCAP)-1 with its putative receptor, Latrophilin-1 (ADGRL1), in immortalized cell lines". University of Toronto, M.S. thesis, p. 28 (2016).
7. Farahzad, Ali, "Role of a SET/Smad7 Interaction in Skeletal Myogenesis". York University, M.S. thesis, p. 29 (September 2016).
8. Bignoux, Monique, "*In vivo* and *in vitro* analysis on the effect of LRP/LR on Alzheimer's disease related proteins, TERT expression and telomerase activity in Alzheimer's disease models". University of the Witwatersrand, M.Sc. thesis, p. 71 (2018).
9. Cuttler, Katelyn, "Investigating the Effect of LRP/LR and Telomerase on Tauopathy In Alzheimer's Disease Cell Culture Models". University of the Witwatersrand, M.Sc. thesis, p. 29 (2019).
10. Abdollahzadeh, Elnaz, "Regulation of BMI1-RING1B interaction and its implication in histone modification". California State Polytechnic, Pomona, M.S. thesis, p. 21 (2020).
11. Scharadin, Tiffany M., "TIG3: A multifunctional regulator of cell proliferation and survival". University of Maryland Baltimore, Ph.D. dissertation, p. 92 (2012).
12. Burkard, Christine, "Host Factors Involved in the Entry of Coronaviruses into Mammalian Cells". Universiteit Utrecht, Ph.D. dissertation, p. 74 (2015).
13. Thienel, Constanze, "Sirtuin 1 and Angiotensin II as Inflammatory Modulators in the Development of Diabetes". Universität Basel, Ph.D. dissertation, pp. 53, 60 (2015).
14. Schihada, Hannes, "Novel optical methods to monitor G-protein-coupled receptor activation in microtiter plates" ("Neue optische Methoden zur Messung der Aktivierung von G-Protein-gekoppelten Rezeptoren in Mikrotiter-Platten"). Julius-Maximilians-Universität Würzburg, Ph.D. dissertation, p. 50 (2018).
15. Zhang, Hongsheng, "Identification and Biochemical Characterization of PGC-1 β - Interacting Proteins". University of Kansas, Ph.D. dissertation, p. 20 (December 2018).
16. Toranzo, Ismael Lamas, "Desarrollo y aplicación de la tecnología CRISPR para el estudio de procesos reproductivos en mamíferos" ("Development and application of CRISPR technology for the study of reproductive processes in mammals"). Universidad Complutense de Madrid, Ph.D. dissertation, pp. 56, 118 (2020).

Notice

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

Technical Assistance

Visit the tech service page at [SigmaAldrich.com/techservice](https://www.sigmaaldrich.com/techservice).

Terms and Conditions of Sale

Warranty, use restrictions, and other conditions of sale may be found at [SigmaAldrich.com/terms](https://www.sigmaaldrich.com/terms).

Contact Information

For the location of the office nearest you, go to [SigmaAldrich.com/offices](https://www.sigmaaldrich.com/offices).

Immunofluorescence Troubleshooting Guide

Problem	Possible Cause	Solution
No staining	FLAG® is not expressed on fusion protein.	Verify expression of FLAG® tag on fusion protein by immunoblotting or other method.
	Antibody concentration is not optimal.	<ul style="list-style-type: none"> Determine optimal working dilution for Cy3™ conjugate by titration. Consider using more antibody if no signal or weak signal is detected.
	Inappropriate filter for fluorescent microscopy.	Use the filter recommended for visualizing Cy3™: <ul style="list-style-type: none"> Excitation maximum for Cy3™: 550 nm Emission maximum for Cy3™: 570 nm
	Incubation time with antibody is not adequate.	Increase incubation time.
	Cultured cells: Intracellular expression of FLAG® fusion protein in cultured cells not accessible by antibody.	Cells need to be permeabilized. <ul style="list-style-type: none"> Consider fixing cells in methanol at –20 °C for 10 minutes, followed by 1 minute in acetone at –20 °C. Alternatively, try fixing cells in 3% paraformaldehyde containing 0.5% TRITON® X-100 for 10 minutes at room temperature.
High Background	Aggregates	Centrifuge antibody conjugate briefly in microcentrifuge at highest speed to remove antibody aggregates.
	Antibody binding to Fc receptors on cell surface.	Incubate sample with 10% irrelevant serum, such as goat serum (Cat. No. G9023), to occupy Fc receptors prior to applying the antibody conjugate.
	Wash steps are not adequate.	Increase number or length of washes.
	Antibody concentration is not optimal.	<ul style="list-style-type: none"> Determine optimal working dilution for the Cy3™ conjugate by titration. Consider using less antibody if background is too high.

The life science business of Merck operates as MilliporeSigma in the U.S. and Canada.

Merck, FLAG, ANTI-FLAG, FLAG-BAP, and Sigma-Aldrich are trademarks of Merck or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.

© 2021-2026 Merck and/or its affiliates. All Rights Reserved.

23332169 Rev 01/26

