

CF® Dye Maleimides

CF® Dye Maleimides are thiol-reactive fluorescent dyes. Maleimides are commonly used to label proteins, peptides or other molecules containing free thiol groups.



Product attributes

| | |
|-----------------------------------|--|
| Chemical reactivity (reacts with) | Thiols |
| Functional group | Maleimide |
| Storage Conditions | Store at -10 to -35 °C, Protect from light |

Product Description

CF® Dye Maleimides are thiol-reactive fluorescent dyes. Maleimides are commonly used to label proteins or small ligands like peptides or oligonucleotide thiophosphates on free thiol groups.

- Thiol-reactive fluorescent dyes.
- Stably label antibodies, proteins, or small ligands like peptides or oligonucleotide thiophosphates on free thiol groups.
- Bright, photostable and water-soluble CF® Dyes are excellent options for fluorescent labeling.

Superior CF® Dyes

CF® Dyes are Biotium's line of next-generation fluorescent dyes that have improved brightness, photostability and water solubility compared to other commercially available fluorescent dyes.

Learn more about [CF® Dyes](#). For more information download the [CF® Dye Brochure](#).

CF® Dye Maleimide

| Conjugation | Ex/Em | Size | Catalog No. | Dye Features |
|-------------------------|------------|-----------|-----------------------|----------------------------------|
| CF@350 | 355/450 nm | 1 umol | 92020 | CF@350 Features |
| CF@405S | 411/431 nm | 1 umol | 92030 | CF@405S Features |
| CF@405M | 416/452 nm | 1 umol | 92021 | CF@405M Features |
| CF@405L | 413/547 nm | 1 umol | 92046 | CF@405L Features |
| CF@430 | 424/497 nm | 1 umol | 92118 | CF@430 Features |
| CF@440 | 433/512 nm | 1 umol | 92124 | CF@440 Features |
| CF@450 | 448/533 nm | 1 umol | 96012 | CF@450 Features |
| CF@488A | 490/516 nm | 1 umol | 92022 | CF@488A Features |
| CF@503R | 503/532 nm | 1 umol | 96079 | CF@503R Features |
| CF@532 | 531/552 nm | 1 umol | 92045 | CF@532 Features |
| CF@543 | 543/563 nm | 1 umol | 92044 | CF@543 Features |
| CF@550R | 551/577 nm | 1 umol | 96074 | CF@550R Features |
| CF@555 | 554/568 nm | 1 umol | 92023 | CF@555 Features |
| CF@568 | 562/584 nm | 1 umol | 92024 | CF@568 Features |
| CF@570 | 568/592 nm | 1 umol | 96015 | CF@570 Features |
| CF@583 | 584/606 nm | 1 umol | 96017 | CF@583 Features |
| CF@583R | 585/609 nm | 1 umol | 96107 | CF@583R Features |
| CF@594 | 593/615 nm | 1 umol | 92025 | CF@594 Features |
| CF@620R | 620/643 nm | 1 umol | 92033 | CF@620R Features |
| CF@633 | 629/650 nm | 1 umol | 92026 | CF@633 Features |
| CF@640R | 642/663 nm | 1 umol | 92034 | CF@640R Features |
| CF@647 | 652/668 nm | 1 umol | 92027 | CF@647 Features |
| CF@660R | 663/682 nm | 1 umol | 92031 | CF@660R Features |
| CF@660C | 667/685 nm | 1 umol | 92028 | CF@660C Features |
| CF@680 | 681/698 nm | 1 umol | 92029 | CF@680 Features |
| CF@680R | 680/701 nm | 1 umol | 92032 | CF@680R Features |
| CF@750 | 755/779 nm | 0.5 umol | 96062 | CF@750 Features |
| CF@770 | 770/797 nm | 1 umol | 96132 | CF@770 Features |
| CF@790 | 783/808 nm | 0.25 umol | 96108 | CF@790 Features |
| CF@800 | 797/817 nm | 0.25 umol | 92128 | CF@800 Features |
| CF@820 | 822/835 nm | 0.25 umol | 96069 | CF@820 Features |

CF® Dye Technical Data

| Dye | Abs/Em (nm) | MW (free acid form) | Sephadex® media ¹ | Amax (nm) | Cf A260/Amax | Cf A280/Amax | ε ² | Optimal DOL (IgG) |
|-------------------------|-------------|---------------------|------------------------------|-----------|--------------|--------------|----------------|-------------------|
| CF@350 | 355/450 | ~496 | G-25 | 347 | 0.13 | 0.14 | 18,000 | 4-6 |
| CF@405S | 411/431 | ~1169 | G-25 | 404 | 0.19 | 0.7 | 33,000 | 3-5 |
| CF@405M | 416/452 | ~503 | G-25 | 408 | 0.24 | 0.13 | 41,000 | 4-6 |
| CF@405L | 413/547 | ~1573 | G-25 | 395 | N/A | 0.5 | 24,000 | 8-12 |

| Dye | Abs/Em (nm) | MW (free acid form) | Sephadex® media ¹ | Amax (nm) | Cf A260/Amax | Cf A280/Amax | ε ² | Optimal DOL (IgG) |
|-------------------------|-------------|---------------------|------------------------------|-----------|--------------|--------------|----------------|-------------------|
| CF@410 | 404/455 | ~242 | G-25 | 416 | 0.15 | 0.2 | 46,000 | 5-7 |
| CF@430 | 424/497 | ~429 | G-25 | 426 | 0.21 | 0.044 | 40,000 | 5-8 |
| CF@440 | 433/512 | ~479 | G-25 | 440 | 0.303 | 0.139 | 40,000 | 5-8 |
| CF@450 | 448/533 | ~689 | G-25 | 450 | 0.205 | 0.2 | 40,000 | 5-8 |
| CF@488A | 490/516 | ~914 | G-25 | 490 | 0.16 | 0.1 | 70,000 | 7-9 |
| CF@503R | 503/532 | ~1100 | G-25 | 503 | 0.21 | 0.09 | 90,000 | 4-10 |
| CF@505 | 505/519 | ~587 | G-25 | 505 | 0.22 | 0.09 | 90,000 | 4-8 |
| CF@510 | 513/537 | ~562 | G-25 | 513 | 0.25 | 0.14 | 90,000 | 3-5 |
| CF@514 | 516/549 | ~1216 | G-25 | 516 | 0.14 | 0.073 | 105,000 | 5-8 |
| CF@532 | 531/552 | ~685 | G-25 | 527 | 0.11 | 0.06 | 96,000 | 4-7 |
| CF@543 | 543/563 | ~887 | G-25 | 541 | 0.305 | 0.095 | 100,000 | 4-7 |
| CF@550R | 551/577 | ~686 | G-25 | 551 | 0.12 | 0.08 | 100,000 | 5-6 |
| CF@555 | 554/568 | ~959 | G-25 | 555 | 0.026 | 0.08 | 150,000 | 4-5, 3-6* |
| CF@568 | 562/584 | ~714 | G-25 | 562 | 0.17 | 0.08 | 100,000 | 5-6 |
| CF@570 | 568/592 | ~2998 | G-25 | 568 | 0.0998 | 0.1 | 150,000 | 5-6 |
| CF@583 | 584/606 | ~3127 | G-75 | 583 | 0.139 | 0.223 | 150,000 | 5-6 |
| CF@583R | 585/609 | ~773 | G-25 | 585 | 0.19 | 0.08 | 100,000 | 5-6 |
| CF@594 | 593/615 | ~729 | G-25 | 593 | 0.24 | 0.08 | 115,000 | 4-7 |
| CF@597R | 597/619 | ~800 | G-25 | 597 | 0.25 | 0.08 | 100,000 | 5-6 |
| CF@620R | 620/643 | ~738 | G-25 | 617 | 0.28 | 0.45 | 115,000 | 5-6 |
| CF@633 | 629/650 | ~821 | G-25 | 630 | 0.25 | 0.48 | 100,000 | 4-7 |
| CF@640R | 642/663 | ~832 | G-50 | 642 | 0.23 | 0.44 | 105,000 | 4-7 |
| CF@647 | 652/668 | ~985 | G-25 | 650 | 0.01 | 0.03 | 240,000 | 4-5, 3-6* |
| CF@660C | 667/685 | ~3024 | G-75 | 667 | 0.05 | 0.08 | 200,000 | 3-6, 2-3* |
| CF@660R | 662/682 | ~888 | G-25 | 663 | 0.2 | 0.51 | 100,000 | 4-7 |
| CF@680 | 681/698 | ~3153 | G-75 | 681 | 0.06 | 0.09 | 210,000 | 3-5, 2-3* |
| CF@680R | 680/701 | ~912 | G-25 | 680 | 0.155 | 0.32 | 140,000 | 5-6 |
| CF@700 | 696/721 | ~2474 | G-75 | 696 | 0.055 | 0.06 | 240,000 | 3-6 |
| CF@710 | 713/735 | ~860 | G-25 | 712 | 0.11 | 0.07 | 115,000 | 5-6 |
| CF@725 | 729/750 | ~890 | G-25 | 729 | 0.11 | 0.07 | 120,000 | 5-6 |
| CF@740 | 742/767 | ~900 | G-25 | 742 | 0.132 | 0.08 | 105,000 | 5-6 |
| CF@750 | 755/779 | ~2921 | G-75 | 755 | 0.01 | 0.03 | 250,000 | 3-5, 2-3* |
| CF@770 | 770/797 | ~3091 | G-75 | 770 | 0.041 | 0.06 | 220,000 | 3-5, 2-3* |
| CF@790 | 783/808 | ~3179 | G-75 | 784 | 0.104 | 0.07 | 210,000 | 3-5 |
| CF@800 | 797/817 | ~3334 | G-75 | 797 | 0.09 | 0.08 | 210,000 | 3-5 |
| CF@820 | 822/835 | ~2711 | G-75 | 822 | 0.0459 | 0.07 | 253,000 | 3-6 |
| CF@850 | 852/870 | ~2787 | G-75 | 852 | N/A | 0.06 | 240,000 | 3-6 |
| CF@870 | 876/896 | ~2773 | G-75 | 877 | N/A | 0.06 | 240,000 | 3-6 |

1. Sephadex® media recommendations are for antibody purification, not nucleic acid.

2. Extinction Coefficient (ε).

*Suitable, but suboptimal DOL.

Sephadex is a registered trademark of Cytiva.

References

1. J Gen Physiol (2014) 143 (1): 105–118. [DOI: 10.1085/jgp.201311053](#)
2. J Biol Chem (2015) 290: 17956-17966. [DOI: 10.1074/jbc.M115.641688](#)
3. Nature Methods (2015) 12:773–779. [DOI: 10.1038/nmeth.3475](#)
4. Biophys J (2018) 115:996-1006. [DOI: 10.1016/j.bpj.2018.07.037](#)

Download a list of [CF® Dye references](#).

This datasheet was generated on June 12, 2026 at 12:40:52 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cf-dye-maleimide/>