

CF™405L Dye

A 405-excitable dye with orange fluorescence emission

Technical Summary

Abs/Em Maxima: 395/545 nm

Extinction coefficient: 24,000

Molecular weight: ~ 1573

Flow cytometry laser line: 405 nm

Microscopy laser line: 405 nm

Direct replacement for: Pacific Orange®

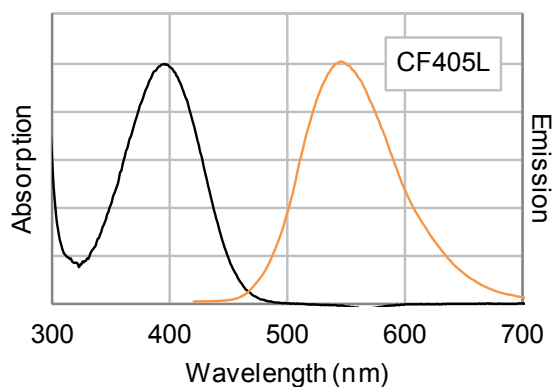


Figure 1. Absorption and emission spectra of CF™405L conjugated to goat anti-mouse IgG in PBS.

CF™405L is a 405-excitable dye with orange fluorescence emission centered at 545 nm (Figure 1). Blue fluorescent CF™405M and orange fluorescent CF™405L can be both excited by the 405 laser and emit at 452 nm and 545 nm, respectively, for multi-color detection in flow cytometry using the 405 nm laser line.

A list of CF™405L-based products is shown in Table 1. A full selection of secondary antibodies, antibody labeling kits, and other bioconjugates including phalloidins, Annexin V and α -bungarotoxin are available for many CF™ dyes. Visit www.biotium.com for details.

CF™405L fluorescent reagents

| Product Name | Size | Cat. No. |
|------------------------------------------------------------------|-------------|-------------|
| CF™405L-Labeled Secondary Antibodies and Other Conjugates | | |
| Goat anti-mouse IgG (H+L), whole antibody, 2 mg/mL | 0.5 mL | 20408-500uL |
| | 50 uL | 20408-50uL |
| | 1 mg | 20408-1mg |
| Goat Anti-Rabbit IgG (H+L) whole antibody, 2 mg/mL | 0.5 mL | 20409-500uL |
| | 50 uL | 20409-50uL |
| | 1 mg | 20409-1mg |
| Streptavidin | 1 mg | 29056 |
| CF™405L Reactive Dyes and Labeling Kits | | |
| Succinimidyl Ester | 1 umol | 92112 |
| Maleimide | 1 umol | 92046 |
| SE Protein Labeling Kit (3 x 1 mg labelings) | 3 labelings | 92228 |
| Mix-n-Stain™ Antibody Labeling Kit (5-20 ug labeling) | 1 labeling | 92303 |
| Mix-n-Stain™ Antibody Labeling Kit (20-50 ug labeling) | 1 labeling | 92304 |
| Mix-n-Stain™ Antibody Labeling Kit (50-100 ug labeling) | 1 labeling | 92305 |

Please visit www.biotium.com for a full listing of CF™405L and other CF™ dye products.

CF is a trademark of Biotium; CF dye technologies are covered by pending US and international patents.

Listed products are for research use only. Not for use in diagnostic or therapeutic procedures. Pacific Orange is registered trademark of Molecular Probes.



Biotium, Inc.

3159 Corporate Place, Hayward, CA 94545

www.biotium.com

US orders: 1-800-304-5357