

CF™620R Dye

A far-red dye excellent for FRET and other applications

Technical Summary

Abs/Em Maxima: 617/639 nm

Extinction coefficient: 115,000

Molecular weight: 738

Direct replacement for: LightCycler® Red 640 (LC640)

Advantages

- Highly water-soluble
- Highly fluorescent
- Extremely photostable

CF™620R dye is a far-red rhodamine-based fluorescent dye. The dye is highly fluorescent and extremely photostable. With its absorption and emission maxima centered at 617 and 639 nm, respectively, the dye may be used as an excellent energy acceptor in fluorescence resonance energy transfer (FRET), or used in multi-color detection applications where the excitation and emission windows can be matched with the spectral profiles of the dye for maximal fluorescence collection. The exceptional water solubility of the dye facilitates bioconjugation in aqueous media and better retain the biological specificity of the conjugates.

Two reactive forms of CF™620R dye are currently available (Table 1). The succinimidyl ester dye is for labeling amino molecules while the maleimide dye is for labeling molecules with a thiol group. For more updated CF620R dye products, please visit Biotium website at www.biotium.com.

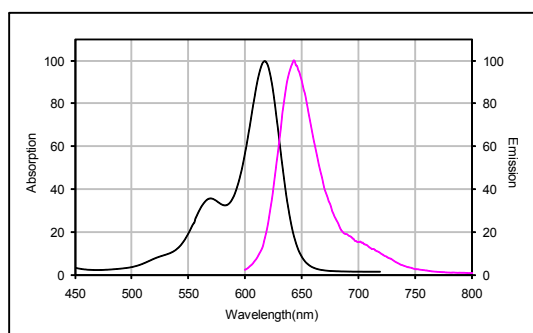


Figure 1. Absorption and emission spectra of CF™620R dye in PBS.

Table 1. CF™620R Product List

Product Name	Size	Cat No.
CF™620R maleimide	1 μmole	92033
CF™620R succinimidyl ester	1 μmole	92106

Listed products are for research use only. Not for use in diagnostic or therapeutic procedures. CF is a trademark of Biotium; CF dye technologies are covered by pending US and international patents. LightCycler is a trademark of Roche.



Biotium, Inc.
3159 Corporate Place, Hayward, CA 94545
www.biotium.com
US orders: 1-800-304-5357