

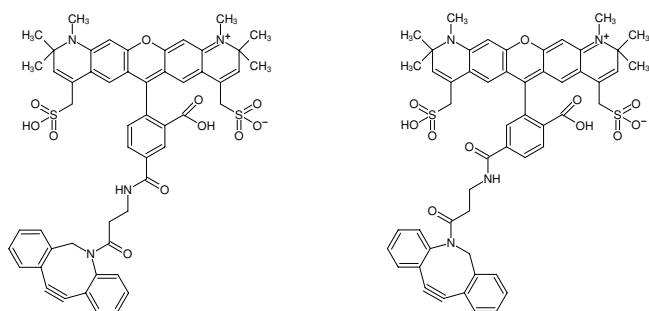


DBC0-AF594

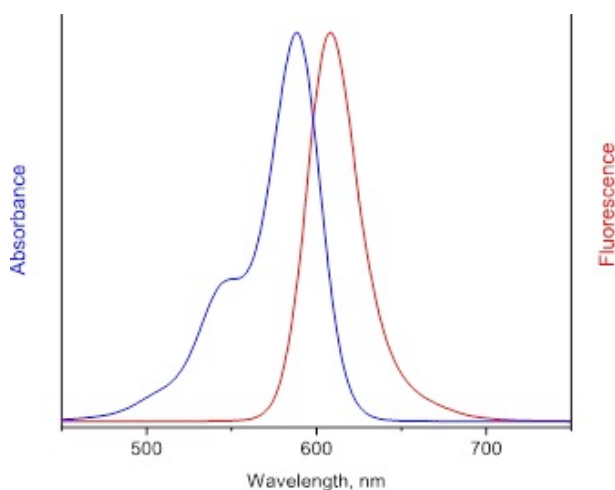
Abs/Em = 590/617 nm

also known as Alexa Fluor® 594-DBC0

| Cat. No. | Amount |
|------------|----------|
| CLK-1298-1 | 1 mg |
| CLK-1298-5 | 5 x 1 mg |



Structural formula of DBC0-AF594



excitation and emission spectrum of AF594

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at -20 °C

Shelf Life: 12 months after date of delivery

Molecular Formula: C₅₃H₄₈N₄O₁₁S₂

Molecular Weight: 981.10 g/mol

Exact Mass: 980.28 g/mol

Purity: ≥ 90 % (HPLC)

Form: solid

Color: blue

Solubility: DMSO, DMF

Spectroscopic Properties: λ_{exc} 590 nm, λ_{em} 617 nm, ε 92.0 L mmol⁻¹ cm⁻¹