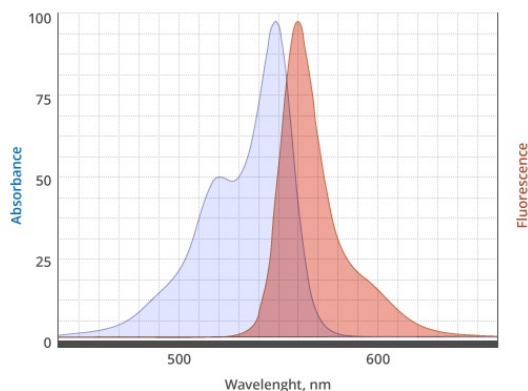


**AF555-Azide**

Abs/Em = 555/572 nm
structural analog to Alexa Fluor® 555

Cat. No.	Amount
CLK-090-1	1 mg
CLK-090-5	5 x 1 mg



excitation and emission spectrum of AF555

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at -20 °C

Shelf Life: 12 months after date of delivery

Molecular Weight: 915.08 g/mol (protonated)

Purity: ≥ 90 % (HPLC)

Form: solid

Solubility: DMSO, DMF

Spectroscopic Properties: λ_{exc} 555 nm, λ_{em} 572 nm,
 ϵ 155.0 L mmol⁻¹ cm⁻¹