



AF594-Picolyl-Azide

Abs/Em = 590/617 nm

also known as Alexa Fluor® 594-Picolyl-Azide

Cat. No.	Amount
CLK-1296-1	1 mg
CLK-1296-5	5 x 1 mg

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: C₄₄H₄₄N₈O₁₁S₂

Molecular Weight: 925.00 g/mol

Exact Mass: 924.26 g/mol

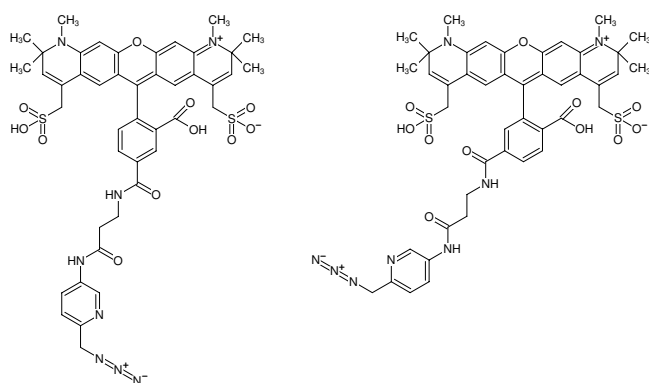
Purity: ≥ 90 % (HPLC)

Form: solid

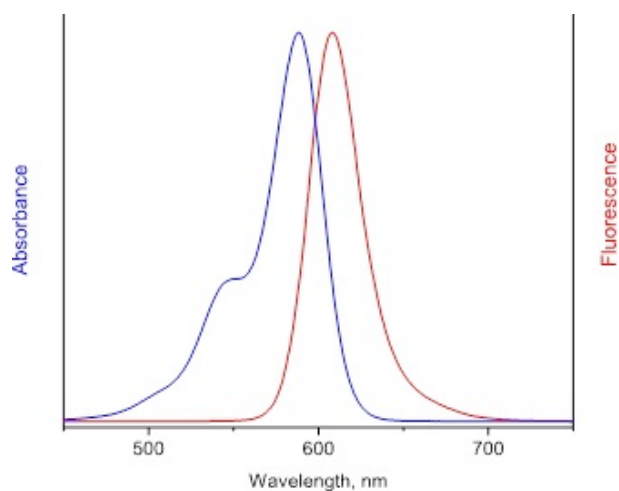
Color: blue

Solubility: DMSO, DMF, MeOH

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



Structural formula of AF594-Picolyl-Azide



excitation and emission spectrum of AF594