



AZDye546-Azide

Abs/Em = 554/570 nm
structural analog to Alexa Fluor® 546

Cat. No.	Amount
CLK-1283A-AZ-1	1 mg
CLK-1283A-AZ-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: C₄₂H₅₃N₇O₁₂S₃

Molecular Weight: 944.10 g/mol

Exact Mass: 943.29 g/mol

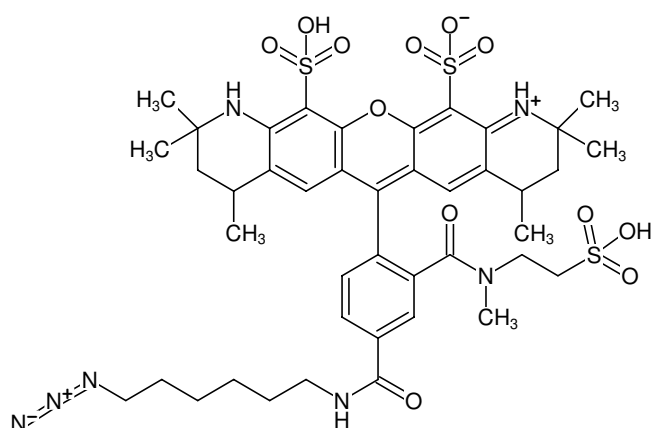
Purity: ≥ 90 % (HPLC)

Form: solid

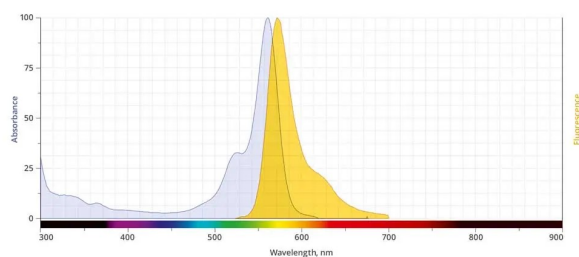
Color: red

Solubility: DMSO, DMF

Spectroscopic Properties: λ_{exc} 554 nm, λ_{em} 570 nm, ε 110.0 L mmol⁻¹ cm⁻¹



Structural formula of AZDye546-Azide



excitation and emission spectrum of AZDye 546