



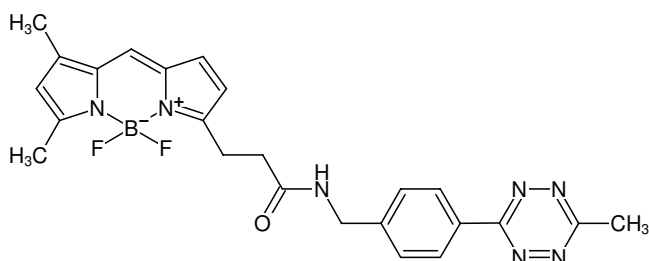
6-Methyl-Tetrazine-BDP-FL

also known as 6-Methyl-Tetrazine-BODIPY® FL

Abs/Em = 503/512 nm

(4-(6-Methyl-1,2,4,5-tetrazin-3-yl)phenyl)methanamine - BDP-FL

Cat. No.	Amount
CLK-037-1	1 mg



Structural formula of 6-Methyl-Tetrazine-BDP-FL

For research use only!

Shipping: shipped on blue ice

Storage Conditions: store at -20 °C

Additional Storage Conditions: store dry and under inert gas

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₂₄H₂₄N₇OBF₂

Molecular Weight: 475.31 g/mol

Exact Mass: 475.21 g/mol

Purity: ≥ 90 % (HPLC)

Form: yellow to orange solid

Solubility: DCM, DMF, DMSO, MeOH

Spectroscopic Properties: λ_{exc} 503 nm, λ_{em} 512 nm, ε 80.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5) (after click reaction)