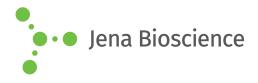
## **DATA SHEET**

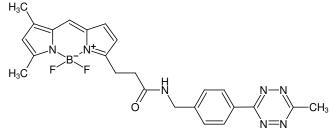




## 6-Methyl-Tetrazine-BDP-FL

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 general laboratory use.

Shipping: shipped on gel packs

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: solid

Color: yellow to orange

Solubility: DCM, DMF, DMSO, MeOH

Spectroscopic Properties:  $\lambda_{exc}$  503 nm,  $\lambda_{em}$  512 nm,  $\epsilon$ 80.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5) (after click reaction)

