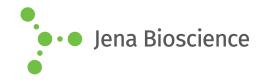
DATA SHEET



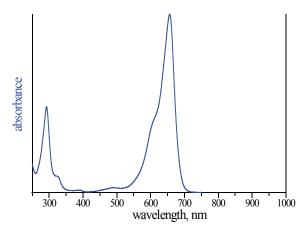


Aminoallyl-dUTP-ATTO-MB2

5-(3-Aminoallyl)-2'-deoxyuridine-5'-triphosphate, labeled with ATTO-MB2, Triethylammonium salt

Cat. No.	Amount
NU-803-MB2	40 μl (1 mM)

Structural formula of Aminoallyl-dUTP-ATTO-MB2



absorption spectrum of ATTO MB2

For general laboratory use.

Shipping: shipped on gel packs **Storage Conditions:** store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient

temperature possible.

Shelf Life: 12 months after date of delivery **Molecular Formula:** C₃₁H₃₉N₆O₁₅P₃S (free acid) **Molecular Weight:** 860.66 g/mol (free acid)

Exact Mass: 860.14 g/mol (free acid)

Purity: ≥ 95 % (HPLC) **Form:** solution in water

Color: blue

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{abs} 668 nm, ϵ 100.0 L mmol⁻¹ cm⁻¹

(Tris-HCl pH 7.5)