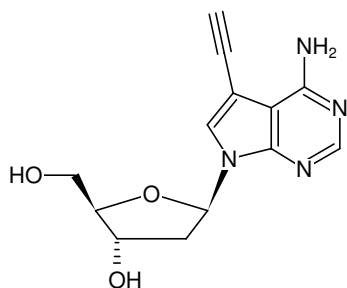


**7-Deaza-7-ethynyl-2'-deoxyadenosine (EdA)**

Cat. No.	Amount
CLK-099	5 mg

Selected References:

Neef *et al.* (2012) Metabolic labeling of DNA by purine analogues in vivo. *Chembiochem.* **13** (12):1750.



Structural formula of 7-Deaza-7-ethynyl-2'-deoxyadenosine (EdA)

For general laboratory use.

Shipping: shipped at ambient temperature

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: C₁₃H₁₄N₄O₃

Molecular Weight: 274.28 g/mol

Exact Mass: 274.11 g/mol

Purity: ≥ 95 % (HPLC)

Form: solid

Color: white to yellow

Spectroscopic Properties: λ_{max} 237/279 nm, ε 11.2/11.1 L mmol⁻¹ cm⁻¹ (Methanol)