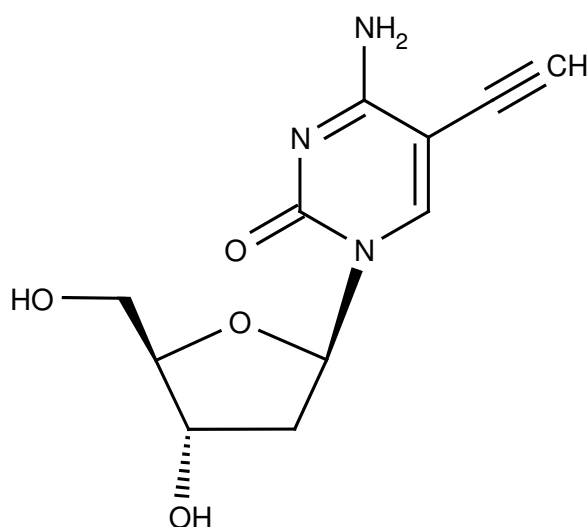


5-Ethynyl-dC (5-EdC) 5-Ethynyl-2'-deoxycytidine

Cat. No.	Amount
CLK-N003-10	10 mg



Molecular Formula: C₁₁H₁₂N₂O₅

Molecular Weight: 252,22 g/mol

Spectroscopic properties: λ_{max} 291 nm; ε 8500 cm⁻¹ M⁻¹ (in Tris-HCl pH 7.5)

Storage conditions: Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

Purity: > 99 % (HPLC)

Appearance: off-white solid

Product Features and Benefits: 5-Ethynyl-2'-deoxycytidine was developed to identify cell proliferation. After incorporation into newly synthesized DNA it can be detected after conjugation to a fluorescent azide via means of click chemistry. It detects DNA at a similar level as 5-Ethynyl-2'-deoxy-uridine (Edu) ¹.

For research use only!

Selected References:

1. Qu *et al.* (2011) 5-Ethynyl-2'-deoxycytidine as a new agent for DNA labeling: detection of proliferating cells. *Anal Biochem* **417(1)**:112.

Guan *et al.* (2011) Intracellular detection of cytosine incorporation in genomic DNA by using 5-ethynyl-2'-deoxycytidine. *Chembiochem* **2(14)**:2184.