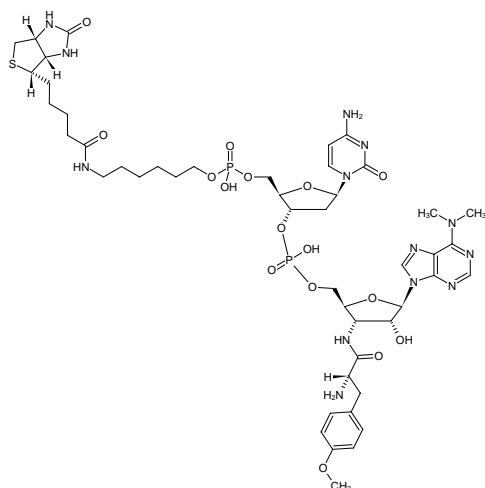




Biotin-dC-puromycin

Cat. No.	Amount
NU-925-BIO-S	100 µl (0,1 mM)
NU-925-BIO-L	5 x 100 µl (0,1 mM)



Structural formula of Biotin-dC-puromycin

For research use only!**Shipping:** shipped on blue ice**Storage Conditions:** store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

Shelf Life: 12 months**Molecular Formula:** C₄₇H₆₉N₁₃O₁₆P₂S (free acid)**Molecular Weight:** 1166.14 g/mol (free acid)**CAS#:** 436083-86-2**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution, pH 7.5 ±0.5**Concentration:** 0.1 mM**pH:** 7.5 ±0.5

Spectroscopic Properties: λ_{max} 260 nm; ε 19.0 L mmol⁻¹ cm⁻¹
(Tris-HCl pH 7.5)

Selected References:

Starck *et al.* (2002) Puromycin oligonucleotides reveal steric restrictions for ribosome entry and multiple modes of translation inhibition. *RNA* **8** (7):890.

Starck *et al.* (2004) A general approach to detect protein expression in vivo using fluorescent puromycin conjugates. *Chem. Biol.* **11** (7):999.

Kawahashi *et al.* (2007) High-throughput fluorescence labelling of full-length cDNA products based on a reconstituted translation system. *J. Biochem.* **141** (1):19.