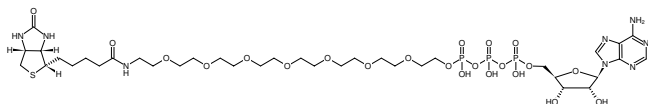


**γ-PEG8-ATP-Biotin**

γ-PEG8-Adenosine-5'-triphosphate-Biotin, Triethylammonium salt

Cat. No.	Amount
NU-926-BIO	50 µl (5 mM)



Structural formula of γ-PEG8-ATP-Biotin

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.

Shelf Life: 12 months after date of delivery**Molecular Formula:** C₃₆H₆₃N₈O₂₂P₃S (free acid)**Molecular Weight:** 1084.91 g/mol (free acid)**Exact Mass:** 1084.30 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 5.0 mM - 5.5 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 259 nm, ε 15.3 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)