

Azide-PEG4-aminoallyl-dUTP

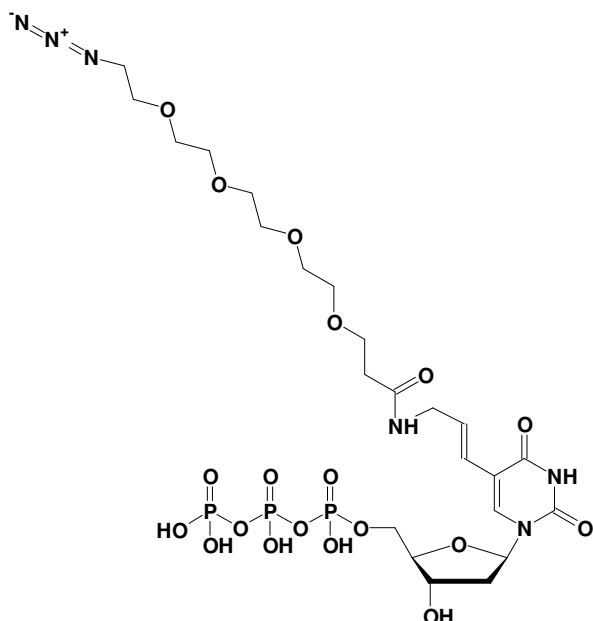
5-(15-Azido-4,7,10,13-tetraoxa-pentadecanoyl-aminoallyl)-2'-deoxy-uridine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-1705S	50 Units
NU-1705L	250 Units

* 1 unit = 1 μ l of a 10 mM solution

Application:

This nucleotide is incorporated into DNA by Taq Polymerase (in-house data).



Molecular Formula: C₂₃H₃₉N₆O₁₉P₃ (free acid)

Molecular Weight: 796.51 (free acid)

Purity: > 95%, clear aqueous solution, pH 7.5

Spectroscopic properties:

λ_{\max} 289 nm; ϵ 7,100 M⁻¹cm⁻¹ 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.

For research use only!