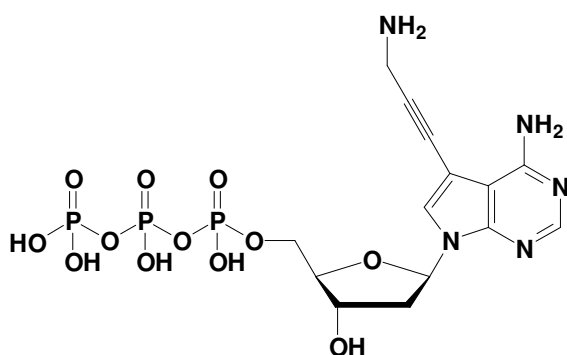


7-Deaza-7-propargylamino-dATP

7-Deaza-7-propargylamino-2'-deoxy-adenosine-5'-triphosphate, sodium salt

Cat. No.	Amount
NU-1611-1	1 mg



Cat. No.: NU-1611

Molecular Formula: $C_{14}H_{20}N_5O_{12}P_3$ (free acid)

Molecular Weight: 543.26 (free acid)

Purity: > 95%

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!