



## 7-Propargylamino-7-deaza-dATP-ATTO-740

7-Deaza-7-propargylamino-2'-deoxyadenosine-5'-triphosphate, labeled with ATTO 740, Triethylammonium salt

Cat. No.	Amount
NU-1611-740	40 µl (1 mM)

**For research use only!**

**Shipping:** shipped on gel packs

**Storage Conditions:** store at -20 °C

**Additional Storage Conditions:** long time storage at pH 7.5

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

**Shelf Life:** 12 months after date of delivery

**Molecular Formula:** C<sub>14</sub>H<sub>20</sub>N<sub>5</sub>O<sub>12</sub>P<sub>3</sub> - ATTO 740 (free acid)

**Molecular Weight:** 992.26 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

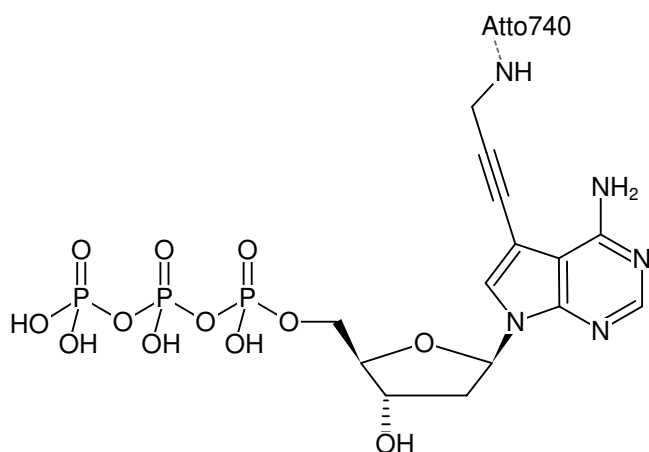
**Form:** solution in water

**Concentration:** 1.0 mM - 1.1 mM

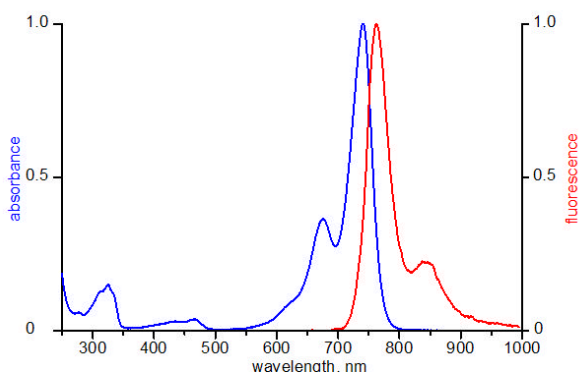
**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 743 nm, λ<sub>em</sub> 763 nm, ε 120.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

**Please note:** The dye decomposes at pH-levels greater 8.0. The recommended pH for a solution of this product is 7.5.



Structural formula of 7-Propargylamino-7-deaza-dATP-ATTO-740



excitation and emission spectrum of ATTO 740