



## 7-Propargylamino-7-deaza-ATP-PEG<sub>5</sub>-AZDye647

structural analog to Alexa Fluor® 647

7-Deaza-7-propargylamino-PEG<sub>5</sub>-adenosine-5'-triphosphate, labeled with AZDye 647, Triethylammonium salt

Cat. No.	Amount
NU-957-PEG5-AZ647	10 µl (5 mM)

**For general laboratory use.**

**Please centrifuge briefly before opening (volume ≤2 ml).**

**Shipping:** shipped on gel packs

**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:** C<sub>64</sub>H<sub>91</sub>N<sub>8</sub>O<sub>32</sub>P<sub>3</sub>S<sub>4</sub> (free acid)

**Molecular Weight:** 1705.62 g/mol (free acid)

**Exact Mass:** 1704.38 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

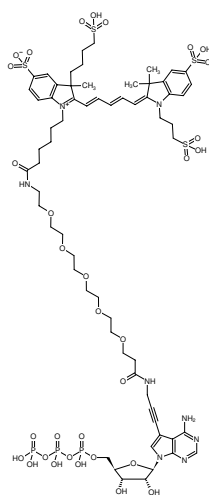
**Form:** solution in water

**Color:** blue

**Concentration:** 5.0 mM - 5.5 mM

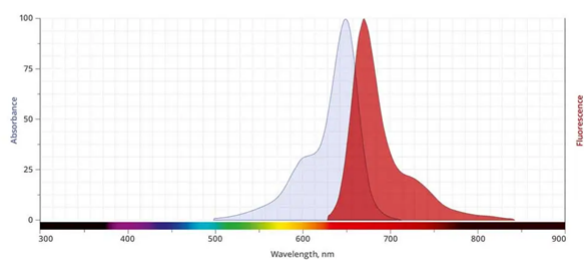
**pH:** 7.5 ±0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 649 nm, λ<sub>em</sub> 671 nm, ε 270.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)



Structural formula of 7-Propargylamino-7-deaza-ATP-PEG<sub>5</sub>-AZDye647

Abs/Em Spectra



excitation and emission spectrum of AZDye 647