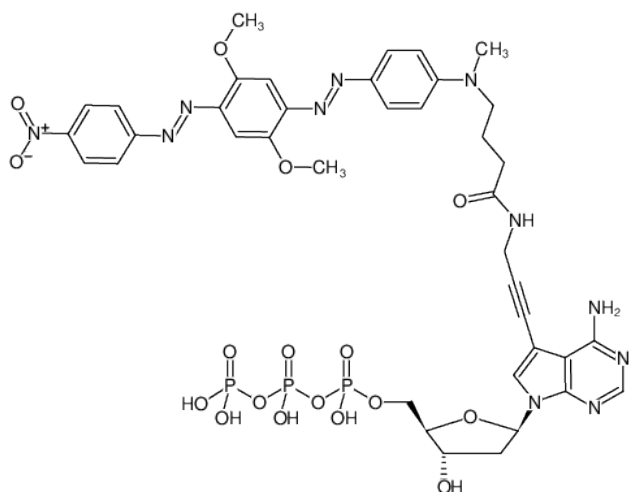




### 7-Propargylamino-7-deaza-dATP-BHQ-2

7-Deaza-7-propargylamino-2'-deoxyadenosine-5'-triphosphate, labeled with BHQ-2™, Triethylammonium salt

Cat. No.	Amount
NU-1611-BHQ2	30 µl (1 mM)



Structural formula of 7-Propargylamino-7-deaza-dATP-BHQ-2

#### For general laboratory use.

**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>39</sub>H<sub>44</sub>N<sub>11</sub>O<sub>17</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 1031.76 g/mol (free acid)

**Exact Mass:** 1031.21 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

**Color:** dark blue

**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ±0.5

**Spectroscopic Properties:** λ<sub>max</sub> 579 nm, ε 38.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)