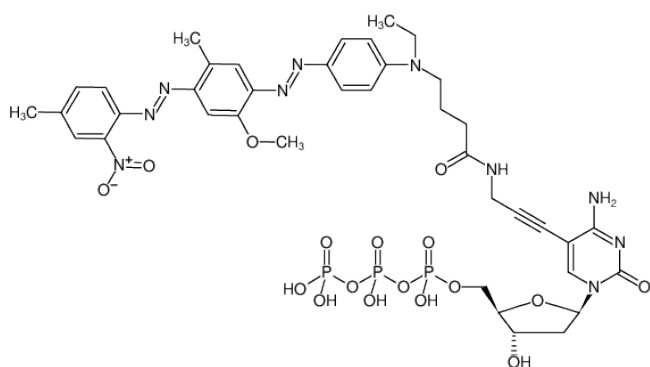




### 5-Propargylamino-dCTP-BHQ-1

5-Propargylamino-2'-deoxycytidine-5'-triphosphate, labeled with BHQ-1™, Triethylammonium salt

Cat. No.	Amount
NU-809-BHQ1	30 µl (1 mM)



Structural formula of 5-Propargylamino-dCTP-BHQ-1

**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>47</sub>N<sub>10</sub>O<sub>17</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 1020.78 g/mol (free acid)

**Exact Mass:** 1020.23 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

**Color:** dark purple

**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ±0.5

**Spectroscopic Properties:** λ<sub>max</sub> 534 nm, ε 34.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)