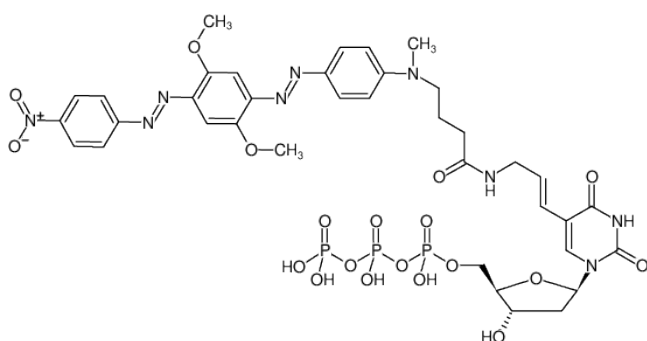




## Aminoallyl-dUTP-BHQ-2

5-(3-Aminoallyl)-2'-deoxyuridine-5'-triphosphate, labeled with BHQ-2™, Triethylammonium salt

Cat. No.	Amount
NU-803-BHQ2	30 µl (1 mM)



Structural formula of Aminoallyl-dUTP-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>37</sub>H<sub>44</sub>N<sub>9</sub>O<sub>19</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 1011.72 g/mol (free acid)

**Exact Mass:** 1011.20 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)