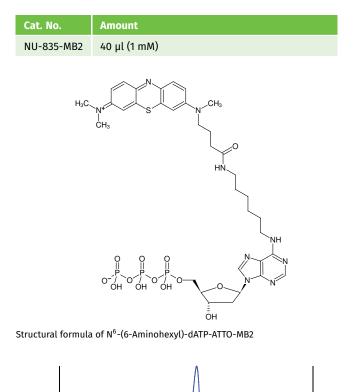




## N<sup>6</sup>-(6-Aminohexyl)-dATP-ATTO-MB2

N<sup>6</sup>-(6-Aminohexyl)-2'-deoxyadenosine-5'-triphosphate, labeled with ATTO-MB2, Triethylammonium salt



## 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>35</sub>H<sub>48</sub>N<sub>9</sub>O<sub>13</sub>P<sub>3</sub>S (free acid)

Molecular Weight: 927.79 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:  $\lambda_{abs}$  668 nm,  $\epsilon$  100.0 L mmol ^1 cm ^1 (Tris-HCl pH 7.5)



400

500

600

700

wavelength, nm

800

900

1000

300

absorbance

