

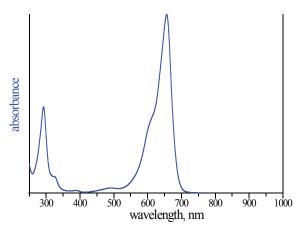


## 8-(6-Aminohexyl)-amino-adenosine-2',5'-bisphosphate-ATTO-MB2

8-(6-Aminohexyl)-amino-adenosine-2',5'-bisphosphate, labeled with ATTO-MB2, Triethylammonium salt

Cat. No.	Amount
NU-812-MB2	40 μl (1 mM)

Structural formula of 8-(6-Aminohexyl)-amino-adenosine-2',5'-bisphosphate-ATTO-MB2



absorption spectrum of ATTO MB2

## 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:  $C_{35}H_{48}N_{10}O_{11}P_2S$  (free acid) Molecular Weight: 878.83 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<sup>-1</sup> cm<sup>-1</sup>

(Tris-HCl pH 7.5)