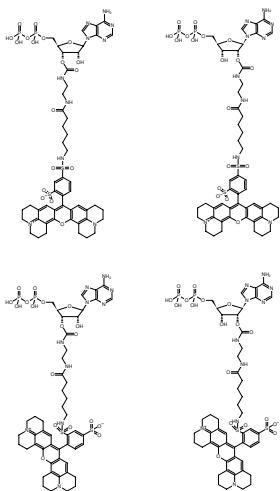


**EDA-ADP-Texas Red**

2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine-5'-diphosphate, labeled with Texas Red, Triethylammonium salt

Cat. No.	Amount
NU-802-TXR	160 µl (1 mM)



Structural formula of EDA-ADP-Texas Red

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: C₅₀H₆₀N₁₀O₁₈P₂S₂ (free acid)

Molecular Weight: 1215.15 g/mol (free acid)

Exact Mass: 1214.30 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: red-violet

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{exc} 588 nm, λ_{em} 609 nm, ε 80.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)