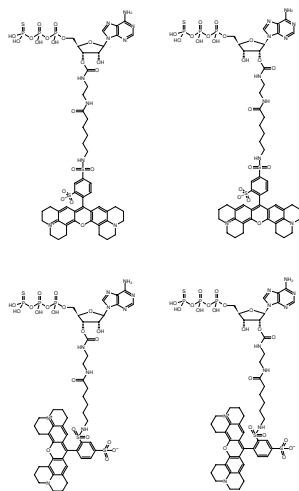




### EDA-ATPyS-Texas Red

2'/3'-O-(2-Aminoethyl-carbamoyl)-adenosine-5'-(γ-thio)-triphosphate, labeled with Texas Red, Triethylammonium salt

Cat. No.	Amount
NU-1609-TXR	50 µl (1 mM)



Structural formula of EDA-ATPyS-Texas Red

#### For general laboratory use.

**Shipping:** shipped on dry ice

**Storage Conditions:** store at -20 °C

**Shelf Life:** 6 months after date of delivery

**Molecular Formula:** C<sub>50</sub>H<sub>61</sub>N<sub>10</sub>O<sub>20</sub>P<sub>3</sub>S<sub>3</sub> (free acid)

**Molecular Weight:** 1311.91 g/mol (free acid)

**Purity:** ≥ 90 % (HPLC)

**Form:** solution in water

**Color:** red-violet

**Concentration:** 1.0 mM - 1.1 mM

**pH:** 7.5 ± 0.5

**Spectroscopic Properties:** λ<sub>exc</sub> 588 nm, λ<sub>em</sub> 609 nm, ε 80.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)