

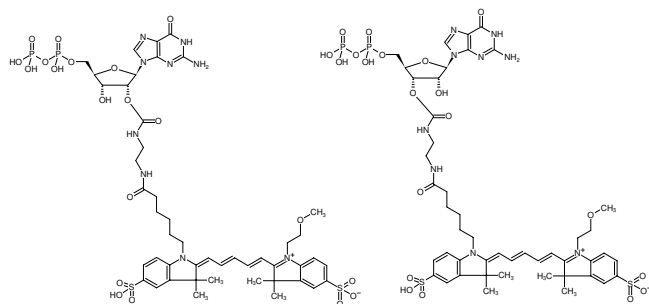
**EDA-GDP-DY-647P1**

2'/3'-O-(2-Aminoethyl-carbamoyl)-guanosine-5'-diphosphate, labeled with DY-647P1, Triethylammonium salt

Cat. No.	Amount
NU-840-647P1	80 µl (1 mM)

Selected References:

Hillig *et al.* (2019) Discovery of potent SOS1 inhibitors that block RAS activation via disruption of the RAS-SOS1 interaction *PNAS* **116** (7):2551.



Structural formula of EDA-GDP-DY-647P1

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: C₄₇H₆₁N₉O₂₀P₂S₂ (free acid)

Molecular Weight: 1198.11 g/mol (free acid)

Exact Mass: 1197.29 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: blue

Concentration: 1.0 mM - 1.1 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{max} 653 nm, ε 250.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)