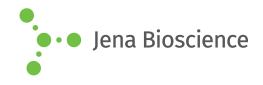
## **DATA SHEET**





## 📕 🔳 γ-[(8-Aminooctyl)-imido]-dUTP

γ-[(8-Aminooctyl)-imido]-2'-deoxyuridine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-818-5	5 mg
NU-818-25	25 mg

H<sub>2</sub>N OH OH OH

Structural formula of  $\gamma$ -[(8-Aminooctyl)-imido]-dUTP

For general laboratory use.

**Shipping:** shipped on gel packs **Storage Conditions:** store at -20 °C

Exact Mass: 594.13 g/mol (free acid)

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

**Shelf Life:** 12 months after date of delivery **Molecular Formula:** C<sub>17</sub>H<sub>33</sub>N<sub>4</sub>O<sub>13</sub>P<sub>3</sub> (free acid) **Molecular Weight:** 594.39 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

Form: solid

Color: white to off-white

**Please note:** This compound contains a phosphoramide linkage which is hydrolyzed at pH <7.0.

## **Selected References:**

Suto *et al.* (1998) Synthesis of y-phosphate-linked nucleoside affinity chromatography resins for protein purification, including ribonucleoside triphosphate reductase. *Nucleosides and Nucleotides* **17 (8)**:1453.