

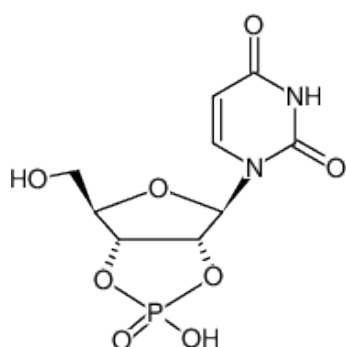


## Uridine-2',3'-cyclic monophosphate

2',3'-cUMP

Uridine-2',3'-cyclic monophosphate, Monosodium salt

Cat. No.	Amount
NU-922-50	50 mg
NU-922-250	5 x 50 mg



Structural formula of Uridine-2',3'-cyclic monophosphate

**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<sub>9</sub>H<sub>11</sub>N<sub>2</sub>O<sub>8</sub>P (free acid)

**Molecular Weight:** 306.17 g/mol (free acid)

**Exact Mass:** 306.03 g/mol (free acid)

**CAS#:** 15718-50-0

**Purity:** ≥ 95 % (HPLC)

**Form:** solid

**Color:** white to off-white

**Spectroscopic Properties:** λ<sub>max</sub> 262 nm, ε 9.8 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

### Selected References:

Marotta *et al.* (2023) Insights into the metabolism, signaling, and physiological effects of 2',3'-cyclic nucleotide monophosphates in bacteria. *Critical Reviews in Biochemistry and Molecular Biology*. **58(2-6)**:118.

Chauhan *et al.* (2022) Binding of 2',3'-Cyclic Nucleotide Monophosphates to Bacterial Ribosomes Inhibits Translation. *ACS Cent. Sci.* **8(11)**:1518.

Duggal *et al.* (2022) Cellular Effects of 2',3'-Cyclic Nucleotide Monophosphates in Gram-Negative Bacteria. *ACS Cent. Sci.* **204(1)**:e00208-21.