



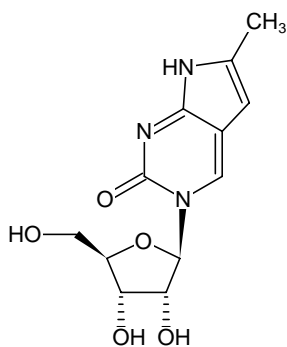
## Pyrrolo-C

6-Methyl-3-(β-D-2-ribofuranosyl)pyrrolo[2,3-d]pyrimidin-2-one

Pyrrolo-C

6-Methyl-3-(beta-D-2-ribofuranosyl)pyrrolo[2,3-d]pyrimidin-2-one

Cat. No.	Amount
N-PYA11092-100	100 mg
N-PYA11092-1G	1 g



Structural formula of Pyrrolo-C

**For research use only!**

**Shipping:** shipped on blue ice

**Storage Conditions:** store at 4 °C

**Shelf Life:** 12 months

**Molecular Formula:** C<sub>12</sub>H<sub>15</sub>N<sub>3</sub>O<sub>5</sub>

**Molecular Weight:** 281.27 g/mol

**CAS#:** 644962-84-5

**Purity:** ≥ 97 %

**Form:** solid

### Selected References:

Berry *et al.* (2004) Pyrrolo-dC and pyrrolo-C: fluorescent analogs of cytidine and 2'-deoxycytidine for the study of oligonucleotides. *Tetrahedron Lett* **45** (11):2457.

Thompson *et al.* (2005) Properties of a New Fluorescent Cytosine Analogue, Pyrrolocytosine. *J. Phys. Chem. B* **109**:6012.