



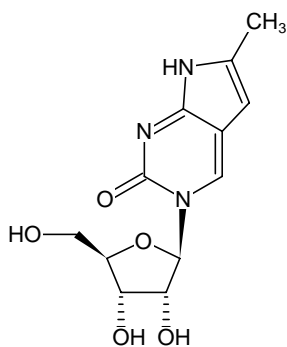
Pyrrolo-C

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 at ambient temperature

Storage Conditions: store at 4 °C

Shelf Life: 24 months after date of delivery

Molecular Formula: C₁₂H₁₅N₃O₅

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.