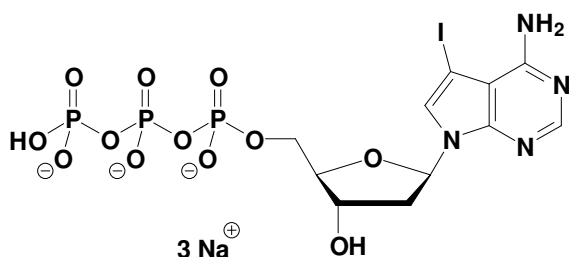


## 7-Deaza-7-I-dATP

7-Deaza-7-Iodo-2'-Deoxy-Adenosine-5'-triphosphate, Sodium salt

| Cat. No. | Amount    |
|----------|-----------|
| NU-1181S | 50 Units  |
| NU-1181L | 250 Units |



**Cat. No.:** NU-1181

**Molecular Formula:** C<sub>11</sub>H<sub>13</sub>N<sub>4</sub>O<sub>12</sub>P<sub>3</sub>I (Anion)

**Molecular Weight:** 613.06 (Anion)

**Purity:** > 95%, clear aqueous solution, pH 7.5

**Storage conditions:**

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

\* 1 unit = 1 µl of a 10 mM solution

**For research use only!**