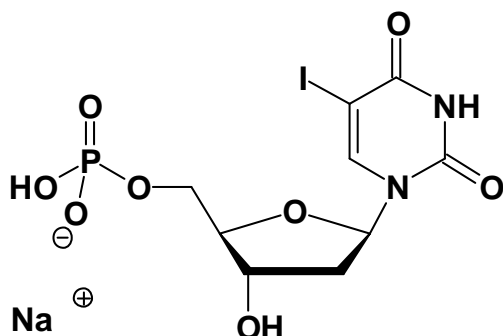


## 5I-dUMP

5-Iodo-2'-deoxy-uridine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-130S	100 Units
NU-130L	500 Units



**Cat. No.:** NU-130

**Molecular Formula:** C<sub>9</sub>H<sub>11</sub>N<sub>2</sub>O<sub>8</sub>PI (Anion)

**Molecular Weight:** 433.07 (Anion)

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

**Spectroscopic properties:**

$\epsilon_{\max} = 7700 \text{ M}^{-1} \text{ cm}^{-1}$  at 287 nm (Tris-HCl 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.

**For research use only!**

1 unit = 1  $\mu\text{l}$  of a 10 mM solution