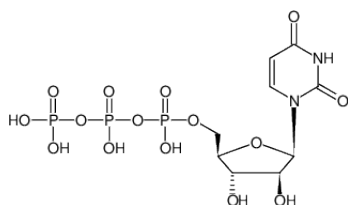



**ara-Uridine-5'-triphosphate (ara-UTP)**

Uridine-arabinofuranoside-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1218S	50 µl (10 mM)
NU-1218L	5 x 50 µl (10 mM)


 Structural formula of *ara*-Uridine-5'-triphosphate (*ara*-UTP)

**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>15</sub>N<sub>2</sub>O<sub>15</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 484.14 g/mol (free acid)

**Exact Mass:** 483.97 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

**Color:** colorless to slightly yellow

**Concentration:** 10 mM - 11 mM

**pH:** 7.5 ±0.5

**Spectroscopic Properties:** λ<sub>max</sub> 262 nm, ε 10.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)