

**Sodium Acetate Solution 3 M pH 5.2**

Cat. No.	Amount
BU-113	100 ml

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at 4 °C

Shelf Life: 12 months

Molecular Formula: CH₃COONa

Molecular Weight: 82.03 g/mol

CAS#: 127-09-3

Form: liquid

Concentration: 3 M

pH: 5.2

Description:

Sodium Acetate Solution has a buffering range from pH 3.6 - 5.6. It is used in the purification and precipitation of nucleic acids, protein crystallization, staining of gels in protein electrophoresis and HPLC.

Quality:

- Molecular biology grade
- sterile filtered, filter pore size 0.22 µm
- DNase / RNase free

