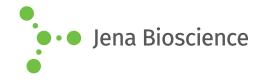
DATA SHEET





■ Potassium Acetate Solution 3 M pH 5.2

Cat. No.	Amount
BU-107	100 ml

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at 4 °C

Shelf Life: 12 months

Molecular Formula: CH₃COOK **Molecular Weight:** 98.14 g/mol

CAS#: 127-08-2 **Form:** liquid

Concentration: 3 M **pH:** 5.2 ±0.1 (22 °C)

Description:

Potassium Acetate Solution has a buffering range from pH 3.8 - 5.8. It is used in the purification and precipitation of nucleic acids. The Buffer is also used to precipitate dodecyl sulfate (DS) and DS-bound proteins allowing the removal of proteins from DNA.

Quality:

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