DATA SHEET





Potassium Chlorid Solution 3 M

Cat. No.	Amount
BU-108-3M	100 ml

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at 4 °C

Shelf Life: 12 months **Molecular Formula:** KCl

Molecular Weight: 74.55 g/mol

CAS#: 7447-40-7 **Form:** liquid

Concentration: 3 M

Description:

Potassium Chloride may be used for the preparation of phosphate buffered saline. It is a component of polymerase chain reaction (PCR) buffer. KCl is also utilized in the extraction, purification and solubilization of proteins.

Quality:

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