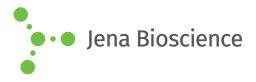
## **DATA SHEET**





## Magnesium Chloride Solution 1 M

Cat. No.	Amount
BU-110-1M	100 ml

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at 4 °C

Shelf Life: 12 months

Molecular Formula: MgCl<sub>2</sub>

Molecular Weight: 95.21 g/mol

CAS#: 7786-30-3

Form: liquid

Concentration: 1 M

## Description:

Magnesium Chloride is used as a source of magnesium ions and for the preparation of competent cells for transformation. It is an essential Co-factor in many enzymes (DNase) and it is included in the polymerase chain reaction (PCR) buffers.

## Quality:

- Molecular biology grade
- sterile filtered, filter pore size 0.22  $\mu m$
- DNase / RNase free

