

**MgCl₂ Solution**

Cat. No.	Amount
PCR-266-25	4 x 1,5 ml (25 mM)

For in vitro use only!

Shipping: shipped at ambient temperature

Storage Conditions: store at 4 °C or -20 °C

Shelf Life: 12 months

Form: liquid

Concentration: 25 mM

Description:

Ultrapure and sterile MgCl₂ solution recommended for individual optimization of PCR assays.

Content:**MgCl₂ (yellow cap)**

25 mM MgCl₂ in PCR-grade water