

**MgCl₂ Solution - 100 mM**MgCl₂ Solution for PCR assay optimization

Cat. No.	Amount
PCR-282S	4 x 1,5 ml (100 mM)
PCR-282-10ML	10 ml (100 mM)
PCR-282-100ML	100 ml (100 mM)

Description:Ultrapure and sterile filtered MgCl₂ solution recommended for individual optimization of PCR assays.**Content:****MgCl₂ Solution**100 mM MgCl₂ in PCR-grade water**For general laboratory use.****Shipping:** shipped at ambient temperature**Storage Conditions:** store at 4 °C or -20 °C**Shelf Life:** 12 months**Form:** liquid**Concentration:** 100 mM