

# 5x siRNA Buffer

This protocol is for 100 mL of 5x siRNA Buffer

## Materials Required:

### Consumables

1. 100 mL sterile bottle or flask
2. 15 mL conical tube

### Chemicals

1. RNase-free water
2. Potassium chloride
3. HEPES (free acid)
4. Magnesium chloride 6H<sub>2</sub>O
5. Potassium hydroxide pellets

### Procedure

1. Prepare 5x siRNA buffer in a sterile 100 mL tube:
  - a. 30 mL RNase-free water
  - b. 2.24 g potassium chloride
  - c. 0.72 g HEPES (free acid)
  - d. 0.02 g magnesium chloride 6H<sub>2</sub>O
2. Prepare fresh 2 M KOH in a separate 15 mL conical tube:
  - a. 1.2 g potassium hydroxide pellets
  - b. 10 mL of RNase-free water
3. Use the 2 M KOH for adjusting the pH of the 5x siRNA buffer from step 1
4. The pH should be between 7.3-7.6
5. Add sterile RNase-free water and mix to a final volume of 100 mL
6. Sterile filter or autoclave the solution. Store the solution at 4 °C for up to 12 months.

GE, imagination at work and GE monogram are trademarks of General Electric Company. Dharmacon is a trademark GE Healthcare companies. All other trademarks are the property of General Electric Company or one of its subsidiaries. ©2014 General Electric Company—All rights reserved. Version published June 2014. GE Healthcare UK Limited, Amersham Place, Little Chalfont, Buckinghamshire, HP7 9NA, UK

