

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 6H₂O
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 6H₂O
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.

If you have any questions

t +44 (0) 1223 976 000 (UK) or +1 800 235 9880 (USA); +1 303 604 9499 (USA)

f + 44 (0)1223 655 581

w horizondiscovery.com/contact-us or dharmacon.horizondiscovery.com/service-and-support

Horizon Discovery, 8100 Cambridge Research Park, Waterbeach, Cambridge, CB25 9TL, United Kingdom

©2018 Horizon Discovery Group Company—All rights reserved. First published April 2018. UK Registered Head Office: Building 8100, Cambridge Research Park, Cambridge, CB25 9TL, United Kingdom.