20X SSC

Saline-Sodium Citrate Salts

Saline-Sodium Citrate (SSC) Concentrate can be used in protocols for Southern blotting, *in situ* hybridization, DNA Microarray or Northern blottings. SSC provides high ionic strength to optimize the rate of probe annealing to the target nucleic acid. SSC is used in hybridization cocktails containing DNA and RNA probes for the analysis of unknown sequences and to locate desired genes and gene fragments. SSC is also used to control stringency for washing steps during *in situ* hybridization. Packet of salts yields 250mL of 20X SSC buffer.

Ordering Information

Order Online »

ENZ-GEN426-0250

250ml

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Formulation Lyophilized.

Reconstitution Dissolve entire content of packet in 250mL of distilled or

deionized water, makes 20X SSC solution.

Last modified: May 29, 2024

