

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.

Citations: 1

[View Online »](#)

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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com