

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.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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