

dsRNA 142 bp

Positive control for dsRNA detection by SCICONS Anti-dsRNA monoclonal J2, K1 und K2 antibodies

SCICONS dsRNA 142 bp is a chemically defined 142 bp long dsRNA sequence intended to be used as a positive control in double-stranded (ds)RNA detection. The sequence does not contain any viral or potentially transforming elements. It is suitable for use as a size standard in dsRNA and RNAi analysis on both non-denaturing polyacrylamide and agarose gels. We suggest to use it as a homogenous, well-characterized positive control for different dsRNA detection methods, e.g. for ELISA and dot-blots.

Ordering Information

[Order Online »](#)

JBS-RNT-SCI-10080100	50µg
----------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Shipping

Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



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