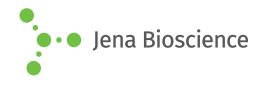
DATA SHEET





■ dsRNA 142 bp

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

Cat. No.	Amount
RNT-SCI-10080100	50 µg

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C to -80 °C upon reconstitution

Additional Storage Conditions: avoid freeze/thaw cycles, store in

aliquots

Shelf Life: 12 months

Form: lyophilised from a 1 mg/ml solution in STE

Solubility: To prepare a 1 μ g/ μ l solution add 50 μ l RNAse-free H₂O.

Description:

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.

Quality control:

Purity: absorbance spectrum and gel electrophresis Activity: Sandwich-ELISA (relative activity to reference antigen)

