

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

DNA for Hybridization Solutions

Catalog #	Contents	Size
9605-5-D	Herring Sperm DNA, Sheared	5 x 1 ml

Description: Highly purified for hybridization (Northern, Southern, probe array) and dot blotting procedures, Trevigen's DNAs are provided at 10 mg/ml with a fragment size range of 200 - 500 bp. The format allows convenient addition to buffers to create a final recommended working concentration of 0.1 mg/ml.

Performance: Quality control testing includes concentration determination, fragment sizing by gel electrophoresis, A_{260}/A_{280} ultraviolet absorbance ratio, and protein content.

Storage Conditions: Store at -20°C in a manual defrost freezer. For long term storage, freeze in working aliquots at -80°C. Avoid repeated freeze-thawing.

References: Meinkoth, J. and Wahl, G. 1984. Hybridization of nucleic acids immobilized on solid supports. *Anal. Biochem.* **138**:267-284 .

Lankford, A.R., Byford, A.M., Ashton, K.J., French, B.A., Lee, J.K., Headrick, J.P. and Matherne, G.P. 2002. Gene expression profile of mouse myocardium with transgenic overexpression of A1 adenosine receptors. *Physiol. Genomics* 11:81-89.

Yeakley, J. M., Fan, J-B., Doucet, D., Luo, L., Wickham, E., Ye, Z., Chee, M.S. and Fu, X-D. 2002. Profiling alternative splicing on fiber-optic arrays. *Nat. Biotechnol.* 20:353-358.

Related Products:

Catalog #	Description	Size
9610-5-D	Salmon Sperm DNA	5 x 1 ml
9600-5-D	Calf Thymus DNA	5 x 1 ml

TREVIGEN®

8405 Helgerman Court, Gaithersburg, MD 20877 USA

Voice: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800

Fax: 301-560-4973 • e-mail: info@trevigen.com • www.trevigen.com

**Herring Sperm DNA
(Sheared)**

Catalog #: 9605-5-D

Storage: -20 °C

TREVIGEN®

1-800-873-8443