

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

DNA for Hybridization Solutions

Catalog #	Contents	Size
9600-5-D	Calf Thymus 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.

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

Kahn, RL., Vadigepalli, R., McDonald, MK., Rogers, RF., Gao, GR. and Schwaber, JS. 2008. Dynamic transcriptomic response to acute hypertension in the nucleus tractus solitarius. *Am J Physiol Regul Integr Comp Physiol* **295**:R15-R27

Related Products:

Catalog #	Description	Size
9610-5-D	Salmon Sperm DNA	5 x 1 ml
9605-5-D	Herring Sperm 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

**Calf Thymus DNA
(Sheared)**

Catalog #: 9600-5-D

Storage: -20 °C

TREVIGEN®

1-800-873-8443