

# TREVIGEN® Product Data

*For Research Use Only. Not For Use In Diagnostic Procedures*

## Salmon Sperm DNA

**Catalog #:** 9610-5-D

**Contents:** Salmon Sperm DNA, Sheared

**Size:** 5 x 1 ml

**Description:** Highly purified for hybridization (Northern, Southern, Chromatin Immunoprecipitation, Microarray) and dot blotting procedures, Trevigen's DNAs are provided at 10 mg/ml with a fragment size range of 80 - 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 .

Ortt, K. and Sinha, S. 2009. Chromatin Immunoprecipitation for Identifying Transcription Factor Targets in Keratinocytes. *Methods Mol Biol.* **585**:159-170.

## Related Products:

Catalog #	Description	Size
9605-5-D	Herring 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](mailto:info@trevigen.com) • [www.trevigen.com](http://www.trevigen.com)

**Salmon Sperm DNA  
(Sheared)**

**Catalog #: 9610-5-D**

**Storage: -20 °C**

**TREVIGEN®**

**1-800-873-8443**