

TREVIGEN[®] Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

Activated DNA

Catalog #: 4667-50-06

Size: 500 μ l

Description: Activated DNA is useful as an activator of Poly(ADP-ribose) polymerase. Activated DNA is provided as optimal for PARP activation.

Source: Purified from herring sperm DNA.

Performance: Quality control testing includes concentration determination, A_{260}/A_{280} ultraviolet absorbance ratio, and qualification in Trevigen's Universal PARP Assay.

Storage Buffer: 10 mM Tris-Cl, pH 8.0, 1 mM EDTA.

Storage Conditions: Store at -20°C.

References:

1. Satoh, M.S. and T. Lindahl. 1994. Role of poly(ADP-ribose) formation in DNA repair. *Nature* **356**:356-358.
2. Lautier, D., J. Lagueux, J. Thibodeau, L. Menard, and G.G. Poirier. 1993. Molecular and biochemical features of poly(ADP-ribose) metabolism. *Mol Cell Biochem* **122**:171-193.

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

Activated DNA
Catalog #: 4667-50-06
Storage: -20 °C

TREVIGEN®
1-800-873-8443