
Nucleic Acid Isolation Enhancer (Glycogen Solution)

(Catalog # R-1001)

Size: 500 µl

Product Description:

Used as a carrier for nucleic acid precipitation in ethanol.

Concentration:

20 µg/µl in RNA-free water.

Storage:

Store at 0-4°C. Stable for 1 year from date of shipment.

Usage:

For research use only. Not for diagnostic or therapeutic use.

References:

Galliard C and Strauss F: Nucleic Acid Res, 18: 378, 1990

Nuclei Acid Isolation Enhancer (Glycogen Solution) is for research use only and is not intended for diagnostic or therapeutic application.