
dNTP Mix

10 mM each x 200 μ l

170-8874

For Research purposes only
Store at -20°C

Storage and Stability

Store the dNTP mix at -20°C in a constant temperature freezer. Avoid repeated freeze/thaw cycles. When stored under these conditions the dNTP mix is stable for one year after ship date.

Contents

The mix contains 10 mM of each dNTP (dATP, dCTP, dGTP, and dTTP). Sufficient for 200 x 50 μ l reactions at a final concentration of 200 μ M (use 1 microliter for a 50 microliter reaction)