

Oligonucleotide stability

Eurogentec performed a 2- year stability study on different types of oligonucleotides and different storage conditions. The results of these studies are summarized in the table below. The criterion used to assess the stability was $\leq 5\%$ variation in purity using high-resolution Ion Exchange HPLC and capillary electrophoresis (CGE).

		Storage ¹	Stability ²
Custom Oligonucleotides	Dried	RT	18 months
	TE buffer (pH 8) or dH ₂ O	-20°C	24 months
Real-Time qPCR probes	Dried	RT	18 months
	TE buffer (pH 8) or dH ₂ O	-20°C	24 months
RNAi Oligonucleotides	Dried	RT	18 months
	RNase-free buffer (pH 7.5)	-20 °C	24 months
Catalogue Oligonucleotides	Dried	RT	18 months
PNA Fish Probes / Custom PNA	Dried	RT	18 months

¹ Oligonucleotides and especially dye-labelled oligonucleotides should be protected from light for optimal stability. Tolerance -20°C \pm 5°C.

² Please note that depending on sequence and modifications, the stability of the oligos may vary substantially versus the given values, which should therefore be considered as indicative. Although it is generally recommended to avoid freeze-thaw cycles, we did not observe negative effects in qPCR tests after 10 freeze-thaw cycles.



Robert Spodarek
QC Validation Manager

KANEKA EUROGENTEC S.A.

LIEGE Science Park • Rue Bois Saint-Jean, 5 • 4102 SERAING • BELGIUM
 ☎ +32 4 372 74 00 • 📠 +32 4 372 75 00 • ✉ info@eurogentec.com • 🌐 www.eurogentec.com
 RPM Liège • T.V.A.-(BE)-0427.348.346