

Thermo Scientific Nucleic Acid Technologies

Amidites for Fast Deprotection Conditions:

Preparation and Usage Notes

REAGENT PREPARATION:

To prepare a 0.1 M solution, add the appropriate volume of anhydrous acetonitrile [CH_3CN dried over activated 3 Å molecular sieves (8-12 mesh) for a minimum of 24 hours]. Refer to Table 1 below.

Molecular sieves can be activated by heating overnight under high vacuum at 150-200°C. Once activated, sieves should be stored in a sealed jar, preferably in a desiccator containing a drying agent.

Table 1. Recommended dilution volumes for producing [0.1 M] phosphoramidite solutions

Product Code	Phosphoramidite Quantity (grams)	Volume of solvent to be added (mL)/bottle
27-1723-xx	0.50	5.6
27-1723-xx	1.00	11.3
27-1723-xx	2.00	22.6
27-1723-xx	5.00	56.5
27-1725-xx	0.50	6.3
27-1725-xx	1.00	12.5
27-1725-xx	2.00	62.5
27-1725-xx	5.00	25.0
27-1726-xx	0.50	5.3
27-1726-xx	1.00	10.5
27-1726-xx	2.00	21.1
27-1726-xx	5.00	52.5

OLIGONUCLEOTIDE DEPROTECTION CONDITIONS

Rapid or mild deprotection protocols may be employed when using PAC dA phosphoramidites, IBU dC phosphoramidites or iPr-PAC dG phosphoramidite in conjunction with oligonucleotide synthesis supports and "PAC" amidites. Otherwise, standard deprotection conditions should be used.

Rapid Deprotection Conditions

Incubate at 60°C for 20-60 minutes in concentrated ammonia. If the oligonucleotide is greater than 50 bases in length, incubate for 30-60 minutes at 60°C.

Mild Deprotection Conditions

Incubate at room temperature for 16 hours in concentrated ammonia.

Standard Deprotection Conditions

Incubate at 55°C for 16 hours in concentrated ammonia.

REFERENCES

- Schulhof, J. C., Molko, D., Teoule, R., Nucl. Acids Res. **16**, 319 (1988)
- Schulhof, J. C., Molko, D., Teoule, R., Nucl. Acids Res. **15**, 397 (1987)
- Sinha, N. D., et al., Nucl. Acids Res. **12**, 4539 (1984)
- Sinha, N. D., Biernat, J., Koster, H., Tetrahedron Lett. **24**, 5843 (1983)

Literature code: 0007711E01U