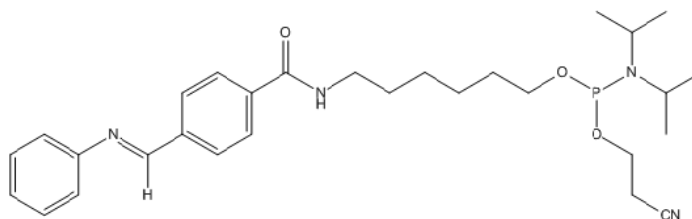


## Product Data Sheet/Certificate of Analysis

### A4FB-Phosphoramidite

Storage: **Store desiccated at or below -20°C**



C<sub>29</sub>H<sub>41</sub>N<sub>4</sub>O<sub>3</sub>P MW 524.63

Catalog Number:	S-1013	Unit Size:	
Lot Number:		Components:	A4FB-Phosphoramidite

Test	Specification	Result
Appearance (color)	Yellow	
Appearance (form)	Oil	
Purity ( NMR)	≥90 % by <sup>32</sup> P NMR	
HPLC analysis after packaging	Conforms to structure	

QC Release Date	Expiration Date:
Released By:	

## Product Description

For generation of 5'-4FB-modified oligonucleotides during solid phase synthesis.

## Comments

**Store desiccated at or below -20°C.** Not for internal or external use in humans.

