

Product Information Sheet

Ordering Information

Product Number: 5260

Product Name: FMOC-Lys(QSY-21)-OH

Unit Size: 50 mg

Storage Conditions:

Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Appearance: Solid

Molecular Weight: 1032.23

Soluble In: DMF

Excitation Wavelength: N/A

Emission Wavelength: N/A

Application Notes

FMOC-Lys(QSY-21)-OH is a building block for in-sequence Lys labeling by QSY-21. QSY-21 is a quencher that can be used to develop a FRET probe with Cy5 or other fluorescent dyes that have the similar fluorescence spectra (to Cy5) such as TF5 or iFluor 647 dyes.