

## FMOC-Lys(QSY-7)-OH

Catalog number: 5261 Unit size: 50 mg

**Product Details** 

Storage Conditions Freeze (<-15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

**Chemical Properties** 

Appearance Solid

Molecular Weight 1032.23

Soluble In DMF

## **Applications**

FMOC-Lys(QSY-7)-OH is a building block for in-sequence Lys labeling by QSY-7. QSY-7 is a quencher that can be used to develop a FRET probe with Cy3, Alexa Fluor 555, iFluor 555, DyLight 555, ATTO 550, ATTO 565 or other fluorescent dyes that have the similar fluorescence spectra (to Cy3) such asTF3 dyes.