

## Product Information Sheet

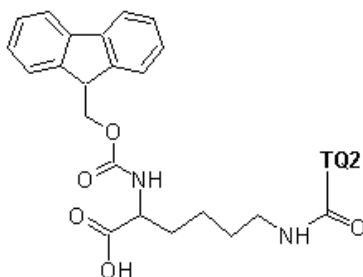
### Ordering Information

Product Number: 5256  
Product Name: FMOC-Lys(TQ2)-OH  
Unit Size: 100 mg  
Storage Conditions: <-15 °C and kept from light and moisture  
Expiration Date: 12 months upon receiving

### Chemical and Spectral Properties

Molecular Weight: 732.89

Chemical Structure:



Soluble in: DMF  
Excitation Wavelength: 515 nm  
Emission Wavelength: NA

### Application Notes

FMOC-Lys(TQ2)-OH is a building block for in-sequence Lys labeling by Tide Quencher™ 2 (TQ2). TQ2 is the best quencher that is used to develop a FRET probe with fluorescein or other fluorescent dyes that have the similar fluorescence spectra (to fluoresceins) such as Bodipy® FL dyes. It can also be used with rhodamine 6G, 6-TET and 6-JOE derivatives.