

TF2-HQKLVFFAK-TQ2 [TF2-His-Gln-Lys-Leu-Val-Phe-Phe-Ala-Lys-TQ2]

Catalog Number: 13206

Unit Size: 1 mg

Product Details

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

Expiration Date 12 months upon receiving

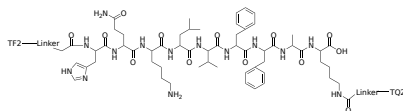
Chemical Properties

Appearance Solid

Molecular Weight N/A

Soluble In DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength 500 nm

Emission Wavelength 522 nm

Applications

TF2-HQKLVFFAK-TQ2 [TF2-His-Gln-Lys-Leu-Val-Phe-Phe-Ala-Lys-TQ2] has much stronger fluorescence upon protease cleavage than the commonly used FRET peptide MCA-HQKLVFFA (K-DNP), a secretase FRET peptide substrate. In addition, the fluorescence of TF2 has much longer wavelength fluorescence than MCA, minimalizing the interference caused by the sample autofluorescence.