



Product Information Sheet

Product Name:	520 MMP FRET Substrate VII
Sequence:	QXL TM 520-Pro-Tyr-Ala-Tyr-Trp-Met-Arg-Lys(5-FAM)-NH ₂ ^{1,2}
Catalog Number:	60574-01
MW:	1965.09
HPLC Purity:	> 95%
Storage:	The product is stable for 1 year at -20°C.

Features and Biological Applications:

This 5-FAM/QXLTM520-based FRET substrate is a sensitive and efficient reagent for assaying MMP activity. It can be cleaved by MMP-7, 12, and 13.

This FRET peptide substrate incorporates QXLTM520, the best quencher available to pair with 5-FAM. When the peptide is intact, the fluorescence of 5-FAM (donor) is quenched by QXLTM520 ("dark" acceptor) through fluorescence resonance energy transfer (FRET). Upon cleavage by MMPs into two separate fragments, the fluorescence of 5-FAM is recovered and can be detected at the emission wavelength of 520±20 nm, with excitation wavelength of 490±20 nm.

Prepare 1 mM DMSO stock solution and dilute in an appropriate assay buffer at a concentration range of 1 to 100 µM. The peptide concentration needs to be optimized depending on your experimental conditions.

References

1. Finch-Arietta, M. et al., *Agents Actions* 39 Spec No, C189 (1993).
2. Netzel-Arnett, S et al. *Anal.Biochem.* 195, 86 (1991).

For *in vitro* research use only