



Product Information Sheet

Product Name:	520 MMP FRET Substrate II
Sequence:	QXL™520 -Pro-Leu-Ala-Leu-Trp-Ala-Arg-Lys(5-FAM)-NH ₂ ¹⁻³
Catalog Number:	60569-01
MW:	1804.1
HPLC Purity:	> 95%
Storage:	The product is stable for 1 year at -20°C.

Features and Biological Applications:

This 5-FAM/QXL™520-based FRET substrate is a sensitive and efficient reagent for assaying MMP activity. It can be cleaved by MMP 1, 7, 8, 12, and 13.

This FRET peptide substrate incorporates QXL™520, the best quencher available to pair with 5-FAM. When the peptide is intact, the fluorescence of 5-FAM (donor) is quenched by QXL™520 (“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. Santala, A. et al. *FEBS Lett.* 461, 153 (1999).
2. Shabani, F. et al. *Free Radic.Res.* 28, 115 (1998).
3. Nagase, H. et al. *J.Biol.Chem.* 269, 20952 (1994).

For *in vitro* research use only