

Product Data Sheet

Product Name: 390 MMP FRET Substrate I

Catalog Number: AS-27076 (1 mg) Lot Number: See label on vial

AS-27077 (5 mg)

Sequence: Mca-Pro-Leu-Gly-Leu-Dap(Dnp)-Ala-Arg-NH2 (3-letter code)

Mca-PLGL-Dap(Dnp)-AR-NH2 (1-letter code)

Molecular Weight: 1094.2

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized yellow powder

Peptide Reconstitution: Using DMSO, reconstitute by adding 80-100 µl to 1 mg MMP FRET

substrate.

Storage: MMP FRET peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at –20°C or lower. Reconstituted peptide can be aliquoted and stored at –20°C or lower.

Description: A fluorogenic substrate that is best cleaved by MMP-7, 8, 9 and 13. The highly fluorescent Mca moiety is efficiently quenched by energy transfer to the Dnp group. The punctuated metallo proteinase (PUMP, EC 3.4.24.23) cleaves the substrate at the Gly-Leu bond with a 190-fold increase in fluorescence (Abs/Em = 325/393 nm). In human matrix metalloproteinases assays, this substrate is about 50 to 100 times more sensitive than the generic MMP substrate, Dnp-PLGLWAr-NH2. Ref: Ref: Knight, CG. et al. *FEBS Lett.* **296**, 263 (1992); Netzel-Arnett, S. et al. *Anal. Biochem.* **195**, 86 (1991).

For Research Use Only