



## Product Data Sheet

---

<b>Product Name:</b>	390 MMP FRET Substrate IV	
<b>Catalog Number:</b>	AS-27109 (5 mg)	Lot Number: See label on vial
<b>Sequence:</b>	Mca-Pro-Cha-Gly-Nva-His-Ala-Dap(Dnp)-NH <sub>2</sub> (3-letter code) Mca-P-Cha-G-Nva-HA-Dap(Dnp)-NH <sub>2</sub> (1-letter code)	
<b>Molecular Weight:</b>	1100.2	
<b>Peptide Purity:</b>	>95%	
<b>Appearance:</b>	Lyophilized yellow powder	

**Peptide Reconstitution:** 390 MMP FRET peptide is freely soluble in water.

**Storage:** 390 MMP FRET peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at  $-20^{\circ}\text{C}$  or lower. Reconstituted peptide can be aliquoted and stored at  $-20^{\circ}\text{C}$  or lower.

**Description:** This substrate is readily hydrolyzed by MMP-8, 9 and collagenase-3 (MMP-13). The hydrolysis is inhibited in a 1:1 stoichiometric fashion by the tissue inhibitors of metalloproteinases, TIMP-1, TIMP-2, and TIMP-3. Abs/Em = 325/393 nm. Ref: Lauer-Fields, JL. and GB. Fields, *Biol. Chem.* **383**, 1095 (2002); Lauer-Fields, JL. et al. *Biochem.* **40**, 5795 (2001); Knauper, V. et al. *J. Biol. Chem.* **271**, 1544 (1996).

### Related Products:

<b>Name</b>	<b>Cat #</b>	<b>Size</b>
390 MMP FRET Substrate I	AS-27076	1 mg
Mca-PLGL-Dpa-AR-NH <sub>2</sub>	AS-27077	5 mg
390 MMP FRET Substrate II	AS-27079	5 mg
Mca-PLGL-Dpa-AR		
390 MMP FRET Substrate V	AS-62027	1 mg
Mca-PLGLEEA-Dpa-NH <sub>2</sub>		

*For Research Use Only*