

Product Data Sheet

Product Name: NFF-3

Catalog Number: AS-27114 (1 mg) Lot Number: See label on vial

Sequence: Mca-Arg-Pro-Lys-Pro-Val-Glu-Nva-Trp-Arg-Lys(Dnp)-NH2 (3-letter code)

Mca-RPKPVE-Nva-WR-K(Dnp)-NH2 (1-letter code)

Molecular Weight: 1675.9

Peptide Purity: >95%

Appearance: Lyophilized yellow powder

Peptide Reconstitution: NFF-3 peptide is freely soluble in water.

Storage: NFF-3 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: NFF-3 is hydrolyzed rapidly by MMP-3 (kcat/Km = 218,000 s-1 M-1) and very slowly by MMP-9 (kcat/Km = 10,100 s-1 M-1), but no significant hydrolysis by MMP-1 and MMP-2. It is one of the few synthetic substrates selectively hydrolyzed by certain members of the MMP family and thus has important applications for the differentiation of MMP-3 activity from that of other MMPs. Abs/Em = 325/393 nm. Ref: Bremer, C. et al. *Acad. Radiol. 9 Suppl* **2**, S314 (2002); Nagase, H. et al. *J. Biol. Chem.* **269**, 20952 (1994).

Published Citations:

Kadonosono, T. et al. Appl. Microbio. Biotech. 75, 1285 (2007).

For Research Use Only