

Beta-Ala-R110

Catalog number: 13203

Unit size: 5 mg

Product Details

Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid
Molecular Weight	545.42
Soluble In	DMSO

Spectral Properties

Excitation Wavelength	500 nm
Emission Wavelength	522 nm

Applications

Beta-Alanine-R110 is a fluorogenic substrate that might be used for detection pancreatic elastase, arylamidase and beta-alanine aminopeptidase activities. It may also be used for detecting *Pseudomonas aeruginosa* that have beta-alanine aminopeptidase activity. Beta-alanine aminopeptidase activity was detected in *Serratia liquefaciens*, *Pseudomonas aeruginosa*, *Pseudomonas fluorescens*, *Burkholderia cepacia*, *Pseudomonas mendocina*., and *Ochrobactrum anthropi*. Therefore, beta-Alanine-R110 may be used for detecting these species in food and other samples. Compared to beta-Alanine-AMC (#13202), beta-Alanine-R110 is more sensitive. Its enzymatic product can be readily detected with the common FITC channel or FITC filter set.