



Product Data Sheet

Product Name: (Z-Ala-Ala-Ala-Ala)2Rh110

Catalog Number: AS-60321-5 (5 mg) **Lot Number:** See label on vial

Sequence: [Rhodamine 110, bis-(CBZ-L-alanyl-L-alanyl-L-alanyl-L-alanine amide)]
(3-letter code)

Molecular Weight: 1170.3

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: (Z-Ala-Ala-Ala-Ala)2Rh110 peptide is freely soluble in DMSO.

Storage: (Z-Ala-Ala-Ala-Ala)2Rh110 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: (Z-Ala-Ala-Ala-Ala)2Rh110 is a very sensitive fluorogenic elastase substrate. The colorless and nonfluorescent (Z-Ala-Ala-Ala-Ala)2Rh110 is hydrolyzed to highly fluorescent rhodamine 110, which exhibits excellent spectral properties that match the optimal detection window of most fluorescence instruments. Abs/Em = 497/520 nm. Ref: Johnson AF, et al. *Anal Chem* **65**, 2352 (1993).

Related Products:

Name	Cat #	Size
(Z-AR)2Rh110•2HCl [Rhodamine 110, bis-(CBZ-L-alanyl-L-arginine amide)]	AS-60322-5	5 mg

For Research Use Only