



Product Data Sheet

Product Name:	Caloxin 2A1	
Catalog Number:	AS-62604 (1mg)	Lot Number: See label on vial
Sequence:	H-Val-Ser-Asn-Ser-Asn-Trp-Pro-Ser-Phe-Pro-Ser-Ser-Gly-Gly-Gly-OH (3-letter code) VSNSNWPSFPSSGGG (1-letter code)	
Molecular Weight:	1479.5	
Peptide Purity:	>95%	
Appearance:	Lyophilized off-white powder	
Peptide Reconstitution:	Caloxin 2A1 is freely soluble in water.	

Storage: Caloxin 2A1 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: Caloxins are extracellular plasma membrane (PM) Ca^{2+} pump inhibitors. Caloxin 2A1 is a PM Ca^{2+} pump inhibitor selectively binding to an extracellular domain. It inhibits Ca^{2+} - Mg^{2+} -ATPase in human erythrocyte leaky ghosts. Caloxin 2A1 is active at an extracellular site, the peptide can simply be added exogenously to inhibit the plasma membrane calcium ATPase (PMCA). Ref: Chaudhary, J. et al. *Am. J. Physiol. Cell. Physiol.* **280**, C1027 (2001); Chaudhary, J. et al. *Am. J. Physiol. Cell. Physiol.* **290**, C1341 (2006); Holmes, M. et al. *Cell Calcium*. **33**, 241 (2003); De Luisi, A. and A. Hofer *J. Cell Sc.* **116**, 1527 (2003).

Related Products:

Name	Cat #	Size
Caloxin 3A1 WSSTSSVSAPLEFGGGGSAK	AS-62606	1 mg

For Research Use Only