



Product Data Sheet

Product Name:	KALA	
Catalog Number:	AS-65459 (1 mg)	Lot Number: See label on vial
Sequence:	Trp-Glu-Ala-Lys-Leu-Ala-Lys-Ala-Leu-Ala-Lys-Ala-Leu-Ala-Lys-His-Leu-Ala-Lys-Ala-Leu-Ala-Lys-Ala-Leu-Lys-Ala-Cys-Glu-Ala-OH WEAKLAKALAKALAKHLAKALAKALKACEA	
Molecular Weight:	3131.8	
Peptide Purity:	Peak Area by HPLC \geq 95%	
Storage:	This 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:	KALA is a cationic amphipathic cell-penetrating peptide (CPP). It assumes an α -helix conformation when the pH is 7.5. KALA binds oligonucleotides and disrupts cell membrane; therefore, it can be used as a DNA transfection reagent.	
References:	<ol style="list-style-type: none">1. Nasrollahi, SA. et al. <i>Chem Biol Drug Des</i> 08, 639-646 (2012).2. Wyman, TB. et al. <i>Biochem</i> 36 (10), 3008-3017 (1997).	

For Research Use Only