



Product Data Sheet

Product Name:	Tat-Beclin-1, scrambled	
Catalog Number:	AS-65468 (1 mg)	Lot Number: See label on vial
Sequence:	Tyr-Gly-Arg-Lys-Lys-Arg-Arg-Gln-Arg-Arg-Arg-Gly-Gly-Val-Gly-Asn-His-Phe-Phe-Ile-Asn-His-Ser-Thr-Thr-Gly-Phe-Ala-Thr-Gln-Trp-OH YGRKKRRQRRRGGVGNDDFFINHETTGFATEW	
Molecular Weight:	3741.1	
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:	Beclin-1 peptide is the HIV-1 Nef binding portion of full-length human Beclin-1 protein (amino acids 267-299). Beclin-1 protein is an autophagy inducing agent that may trigger cellular adaptation, survival, or cell death. When conjugated to the cell-permeable peptide, it can successfully enter cells and induce autophagy. Tat-Beclin-1 (scrambled) will not induce autophagy and can be used as a negative control.	
Related Products:	Tat-Beclin-1, Cat# AS-65467	
References:	<ol style="list-style-type: none">1. Shoji-Kawata, S. et al. <i>Nature</i> 494, 201-209 (2013).2. Marquez, RT. et al. <i>Am J Cancer Res</i> 2 (2), 214–221 (2012).	

For Research Use Only