



## Product Data Sheet

---

<b>Product Name:</b>	Tat-Beclin-1	
<b>Catalog Number:</b>	AS-65467 (1 mg)	Lot Number: See label on vial
<b>Sequence:</b>	Tyr-Gly-Arg-Lys-Lys-Arg-Arg-Gln-Arg-Arg-Arg-Gly-Gly-Thr-Asn-Val-Phe-Asn-Ala-Thr-Phe-Glu-Ile-Trp-His-Asp-Gly-Glu-Phe-Gly-Thr-OH  YGRKKRRQRRRGGTNVFNATFEIWHDGEFGT	
<b>Molecular Weight:</b>	3741.1	
<b>Peptide Purity:</b>	Peak Area by HPLC ≥95%	
<b>Storage:</b>	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.	
<b>Description:</b>	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.	
<b>Related Products:</b>	Tat-Beclin-1, scrambled, Cat# AS-65468	
<b>References:</b>	<ol style="list-style-type: none"><li>1. Shoji-Kawata, S. et al. <i>Nature</i> <b>494</b>, 201-209 (2013).</li><li>2. Marquez, RT. et al. <i>Am J Cancer Res</i> <b>2 (2)</b>, 214–221 (2012).</li></ol>	

*For Research Use Only*