



## Product Data Sheet

<b>Product Name:</b>	Biotin-Exendin 4	
<b>Catalog Number:</b>	AS-60279-05 (0.5 mg) AS-60279-1 (1 mg)	Lot Number: See label on vial
<b>Sequence:</b>	Biotin-His-Gly-Glu-Gly-Thr-Phe-Thr-Ser-Asp-Leu-Ser-Lys-Gln-Met-Glu-Glu-Glu-Ala-Val-Arg-Leu-Phe-Ile-Glu-Trp-Leu-Lys-Asn-Gly-Gly-Pro-Ser-Ser-Gly-Ala-Pro-Pro-Pro-Ser-NH <sub>2</sub> (3-letter code) Biotin-HGEGTFTSDLSKQMEEEEAVRLFIEWLKNGGPSSGAPPPS-NH <sub>2</sub> (1-letter code)	
<b>Molecular Weight:</b>	4412.9	
<b>Peptide Purity:</b>	>95%	
<b>Appearance:</b>	Lyophilized white powder	

**Peptide Reconstitution:** Biotin-Exendin 4 peptide is freely soluble in water.

**Storage:** Biotin-Exendin 4 peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at  $-20^{\circ}\text{C}$  or lower. Reconstituted peptide can be aliquoted and stored at  $-20^{\circ}\text{C}$  or lower.

### Related Products:

<b>Name</b>	<b>Cat #</b>	<b>Size</b>
[Cys(HiLyte Fluor™ 647 C2 maleimide)]-Exendin-4	AS-63714	0.05 mg
C(HiLyteFluor 647 C2 maleimide)-HGEGTFTS DLSKQMEEEEAVRLFIEWLKNGGPSSGAPPPS-NH <sub>2</sub>		
Exendin 4	AS-24463	0.5 mg
HGEGTFTSDLSKQMEEEEAVRLFIEWLKNGGPSSGAPPPS-NH <sub>2</sub>	AS-24464	1 mg
Exendin 4, FAM-labeled	AS-60280-05	0.5 mg
FAM-HGEGTFTSDLSKQMEEEEAVRLFIEWLKNGGPSSGAPPPS-NH <sub>2</sub>		

*For Research Use Only*