



## Product Data Sheet

**Product Name:** Exendin (9-39)  
**Catalog Number:** AS-24467 (0.5 mg)                      Lot Number: See label on vial  
AS-24468 (1 mg)  
**Sequence:** H-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-NH2  
(3-letter code)  
DLSKQMEEEEAVRLFIEWLKNGGPSSGAPPPS-NH2 (1-letter code)  
**Molecular Weight:** 3370.8  
**% Peak Area by HPLC:** ≥ 95  
**Appearance:** Lyophilized white powder  
**Peptide Reconstitution:** Exendin (9-39) peptide is freely soluble in H<sub>2</sub>O.

**Storage:** Exendin (9-39) 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:** This peptide is a potent Glucagon-Like Peptide 1 (GLP-1) receptor antagonist. It is a competitive inhibitor of Exendin-3 and Exendin-4. Ref: Fehmman, HC. et al. *Peptides* **15**, 453 (1994).

### Related Products:

Product Name	Catalog #	Size
Exendin 4	AS-24463	0.5 mg
(HGEGTFTSDLSKQMEEEEAVRLFIEWLKNGGPSSGAPPPS-NH2)	AS-24464	1 mg
Exendin 4, FAM-labeled	AS-60280-05	0.5 mg
(FAM-HGEGTFTSDLSKQMEEEEAVRLFIEWLKNGGPSSGAPPPS-NH2)		
Biotin-Exendin 4	AS-60279-05	0.5 mg
(Biotin-HGEGTFTSDLSKQMEEEEAVRLFIEWLKNGGPSSGAPPPS-NH2)	AS-60279-1	1 mg
Exendin (5-39)	AS-63898	1 mg
(TFTSDLSKQMEEEEAVRLFIEWLKNGGPSSGAPPPS-NH2)		
Fluorescein-Trp25-Exendin-4 (FLEX)	AS-63899	0.5 mg
(HGEGTFTSDLSKQMEEEEAVRLFIE-(Trp-S-FAM)-LKNGGPSSGAPPPS-NH2)		
[Cys(HiLyte™ Fluor 647 C2 maleimide)]-Exendin-4	AS-63714	50 ug
C(HiLyteFluor 647 C2 maleimide)-HGEGTFTSDLSKQMEEEEAVRLFIEWLKNGGPSSGAPPPS-NH2		

*For Research Use Only*