

## **Product Data Sheet**

Product Name:	biotin-Neurogranin (48-76), mouse	
Catalog Number:	AS-65577 (1 mg)	Lot Number: See label on vial
Sequence:	biotin-Ser-Gly-Glu-Cys-Gly-Arg-Lys-Gly-Pro-Gly-Pro-Gly-Gly-Pro-Gly-Gly-Ala-Gly-Gly-Ala-Gly-Gly-Gly-Gly-Gly-Gly-Pro-Ser-OH	
	Biotin-SGECGRKGPGPGGPGGAGGARGGAGGGPS	
Molecular Weight:	2548.9	
Peptide Purity:	Peak Area by HPLC ≥95%	
Storage:	This peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at $-20^{\circ}$ C or lower. Reconstituted peptide can be aliquoted and stored at $-20^{\circ}$ C or lower.	
Description:	Neurogranin is a post synaptic protein that is cytoplasm. It modulates calcium-mediated si critical role in synaptic plasticity. Recently Ne identified as a major neurogranin peptide fra was found that patients with Alzheimer's Dis Neurogranin (48-76) in cerebrospinal fluid.	ignaling pathways and plays eurogranin (48-76) was gment in human brain and it
Related Products:	Neurogranin (48-76), mouse, Cat# AS-65578 Neurogranin (48-76), human, Cat# AS-65578 Biotin-Neurogranin (48-76), human, Cat# AS	5
References:	1. De Vos A., et.al., Alzh and Dement 4, 1-	

- 2. Kvartsberg H., et.al., Alzh Res & Ther 7(40), 1-9 (2015)
- 3. Tarawneh R., et.al., JAMA Neur, E1-E11 (2016)

For Research Use Only