

Product Data Sheet

Product Name: Neurogranin (48-76), human

Catalog Number: AS-65575 (1 mg) Lot Number: See label on vial

Sequence: Ser-Gly-Gly-Arg-Lys-Gly-Pro-Gly-Pro-Gly-Gly-Pro-Gly-Ala-Gly-

Val-Ala-Arg-Gly-Gly-Ala-Gly-Gly-Gly-Pro-Ser-OH

SGERGRKGPGPGGPGGAGVARGGAGGGPS

Molecular Weight: 2417.8

Peptide Purity: Peak Area by HPLC ≥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: Neurogranin is a post synaptic protein that is expressed in dendritic

cytoplasm. It modulates calcium-mediated signaling pathways and plays critical role in synaptic plasticity. Recently Neurogranin (48-76) was identified as a major neurogranin peptide fragment in human brain and it was found that patients with Alzheimer's Disease have elevated levels of

Neurogranin (48-76) in cerebrospinal fluid.

Related Products: Neurogranin (48-76), mouse, Cat# AS-65578

Biotin-Neurogranin (48-76), mouse, Cat# AS-65577 Biotin-Neurogranin (48-76), human, Cat# AS-65576

References:

1. De Vos A., et.al., Alzh and Dement 4, 1-9 (2015)

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