

GAP43 Protein

Full-length recombinant human protein expressed in E. coli cells

Catalog # G15-30H

Lot # B2238-8

Product Description

Full-length recombinant human GAP43 was expressed in E. coli cells using an N-terminal His tag. The gene accession number is [NM_002045](#).

Gene Aliases

B-50; PP46

Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, 25% glycerol.

Storage and Stability

Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

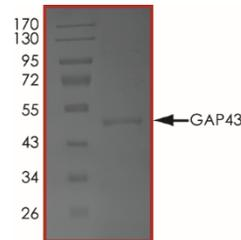
Scientific Background

Growth-associated protein 43 (GAP-43, also known as neuromodulin and B-50) mediates axonal growth, branching, and pathfinding during development. Plays a pivotal role both during development and in axonal remodeling in the adult brain. Increased expression observed in pathological or traumatic lesions induces neuronal rewiring.

References

1. Grasselli G, et al: Structural plasticity of climbing fibers and the growth-associated protein GAP-43. *Front Neural Circuits*. 2013 Feb 21;7:25.
2. Oestreicher A. B., et al: (1997). B-50, the growth associated protein-43: modulation of cell morphology and communication in the nervous system. *Prog. Neurobiol.* 53, 627-686.
3. Buffo A., et al: (2003). Extrinsic regulation of injury/growth-related gene expression in the inferior olive of the adult rat. *Eur. J. Neurosci.* 18, 2146-2158

Purity



The purity of GAP43 was determined to be **>95%** by densitometry. Approx. MW **48 kDa**.

GAP43 Protein

Full-length recombinant human protein expressed in E. coli cells

Catalog #	G15-30H
Lot #	B2238-8
Purity	>95%
Concentration	0.1 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.