

NPR1 Protein

Recombinant human protein expressed in Sf9 cells

Catalog # N22-35G

Lot # M2756-7

Product Description

Recombinant human NPR1 (495-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. This gene accession number is [BC063304](#).

Gene Aliases

ANPa, GUC2A, GUCY2A

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, and 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.

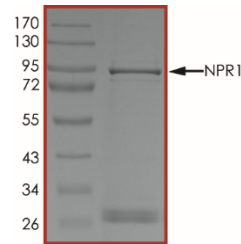
Scientific Background

NPR1 or natriuretic peptide receptor 1 is a membrane-bound guanylate cyclase that serves as the receptor for both atrial and brain natriuretic peptides (1). NPR1 expression influences blood pressures of mice in a dose-dependent manner. ANP and NPR1 binding induces an increase in intracellular cGMP which initiates natriuresis, diuresis, and vasodilation, all of which contribute to lowering blood pressure. Low Npr1 expression leads to a salt-sensitive increase in BP, whereas high Npr1 expression lowers BP and protects against high dietary salt (2).

References

- Garbers, D. L.et.al: Guanylyl cyclase receptors. J. Biol. Chem. 269: 30741-30744, 1994.
- Oliver, P. M.et.al: Natriuretic peptide receptor 1 expression influences blood pressures of mice in a dose-dependent manner. Proc. Nat. Acad. Sci. 95: 2547-2551, 1998.

Purity



The purity of NPR1 protein was determined to be **>75%** by densitometry. Approx. MW **88 kDa**.

NPR1 Protein

Recombinant human protein expressed in Sf9 cells

Catalog #	N22-35G
Lot #	M2756-7
Purity	>75%
Concentration	0.05 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.