

NPR2 Protein

Recombinant human protein expressed in Sf9 cells

Catalog # N23-35G

Lot # M2768-8

Product Description

Recombinant human NPR2 (479-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. This gene accession number is [NM_003995](#).

Gene Aliases

AMDM; ANPB; ANPRB; GUC2B; GUCY2B; NPRB; NPRBI

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.

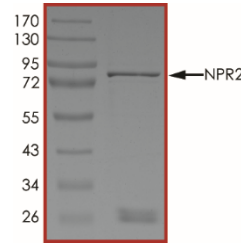
Scientific Background

NPR2 encodes natriuretic peptide receptor B, one of two integral membrane receptors for natriuretic peptides which produce cytoplasmic cyclic GMP from GTP on binding its extracellular ligand, C-type natriuretic peptide (1). The granulosa cell ligand NPPC and its receptor NPR2 in cumulus cells prevent precocious meiotic maturation which is critical for maturation and ovulation synchrony and for normal female fertility. NPR2 is critical for the development of both bone and female reproductive organs (2).

References

1. Lincoln, T. M. et al: Intracellular cyclic GMP receptor proteins. FASEB J. 7: 328-338, 1993.
2. Tamura, N. et al: Critical roles of the guanylyl cyclase B receptor in endochondral ossification and development of female reproductive organs. Proc. Nat. Acad. Sci. 101: 17300-17305, 2004.

Purity



The purity of NPR2 protein was determined to be **>80%** by densitometry. Approx. MW **86 kDa**.

NPR2 Protein

Recombinant human protein expressed in Sf9 cells

Catalog #	N23-35G
Lot #	M2768-8
Purity	>80%
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.