

NF1-333 (RasGAP)

human, recombinant, *E. coli*

Cat. No.	Amount
PR-223	50 µg

For *in vitro* use only
Quality guaranteed for 12 months
Store at -80°C

Avoid freeze / thaw cycles

Form

Liquid. Supplied in 50 mM Tris-HCl pH 7.5, 5 mM MgCl₂ and 3 mM DTE.

Molecular Weight

36.6 kDa (333 amino acids)

Purity

> 95% by SDS-PAGE

Description

The NF1 (Neurofibromin) gene product acts as a GTPase activating factor (GAP) on Ras. Inactivating mutations in NF1 lead to neurofibromatosis type 1.

NF1- 333 contains residues 1198-1530 of human NF1 comprising the functional GAP-related domain that is able to stimulate GTP-hydrolysis on wild type Ras.

Selected References:

Scheffzek *et al.* (1998) Structural analysis of the GAP-related domain from neurofibromin and its implications. *EMBO J.* **17**:4313.

Xu *et al.* (1990) The neurofibromatosis type 1 gene encodes a protein related to GAP. *Cell* **63**:599.

Associated products available from Jena Bioscience

Non-hydrolyzable GTP-analogs (such as GTPγS, GppCp, GppNHp, NPE-caged-GTP...)

Fluorescent GTPs (such as MANT, ANT, TNP, or GTP labeled with a dye of your choice)

Recombinant G-protein signaling reagents (including GAPs, GEFs, heterotrimeric G-proteins, and many more...)

For further information please visit www.jenabioscience.com.