

Ran

Ras-related nuclear protein, TC4

human, recombinant, *E. coli*

Cat. No.	Amount
PR-211	50 µg

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

Avoid freeze / thaw cycles

Form

Liquid. Supplied in 20 mM KH₂PO₄/K₂HPO₄, pH 7.0, 6.5 mM DTE, 150 mM KCl, 5 mM MgCl₂ and 5% glycerol.

Molecular weight

24.0 kDa (216 amino acids).

Purity

> 95% by SDS-PAGE

Description

Ran (Ras-related nuclear protein, TC4) belongs to the superfamily of GTP binding proteins. It acts in diverse cellular processes such as nuclear transport, mitotic spindle assembly, and post-mitotic nuclear assembly. Protein preparation is 90% GDP-loaded, measured by HPLC.

Selected References:

Dasso M (2002) The Ran GTPase: Theme and Variations. *Current Biology* **12**:502.

Renault et al. (2001) Structural Basis for Guanine Nucleotide Exchange on Ran by the regulator of Chromosome Condensation (RCC1). *Cell* **105**:245.

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.