

Ran (E70A)

human, recombinant, *E. coli*

Cat. No.	Amount
PR-215	50 µg

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

Avoid freeze / thaw cycles

Form

Liquid. Supplied as a 1 mg/ml solution in 20 mM $\text{KH}_2\text{PO}_4/\text{K}_2\text{HPO}_4$, pH 7.4, 1 mM DTE and 5 mM MgCl_2 .

Molecular weight

24.0 kDa (216 amino acids).

Purity

> 95% by SDS-PAGE.

Description

Ran (Ras-related nuclear protein, TC4) belongs to the Ras superfamily of GTP binding proteins. It acts in cellular processes such as nuclear transport, mitotic spindle assembly, and post-mitotic nuclear assembly.

The mutation E70A results in inhibition of the guanine exchange reaction mediated by RCC1, the guanine nucleotide exchange factor (GEF) for Ran.

Protein preparation is 85% 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 $\text{GTP}\gamma\text{S}$, 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.