

Ran (Q69L)

human, recombinant, *E. coli*

Cat. No.	Amount
PR-213	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 50 mM $\text{KH}_2\text{PO}_4/\text{K}_2\text{HPO}_4$, pH 7.5, 5 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 superfamily of GTP binding proteins. It acts in diverse cellular processes such as nuclear transport, mitotic spindle assembly, and post-mitotic nuclear assembly. Ran Q69L is defective in GTP hydrolysis. Useful as a negative control for Ran binding and import studies.

Protein preparation is 34% GDP- and 17% GTPloaded, measured by HPLC.

Selected References:

- Dasso (2002) The Ran GTPase: Theme and Variations. *Current Biology* **12**:502.
Palacios *et al.* (1996) Ran/TC4 mutants identify a common requirement for snRNP and protein import into the nucleus. *J. Cell Biol.* **133**:485.
Clarke *et al.* (1995) Regulation of Cdc2/cyclin B activation by Ran, a Ras-related GTPase. *J. Cell Sci.* **108**:1217.

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.