

**Ran - T24N**

Ras-related nuclear protein, TC4
human, recombinant, *E. coli*

Cat. No.	Amount
PR-212	50 µg

For in vitro use only!

Shipping: shipped on dry ice

Storage Conditions: store at -80 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Molecular Weight: 24.0 kDa (216 amino acids)

Accession number: NM_006325

Purity: > 90 % (SDS-PAGE)

Form: liquid (Supplied in 20 mM KH₂PO₄/K₂HPO₄ pH 7.0, 6.5 mM DTE, 150 mM KCl, 5 mM MgCl₂ and 10% glycerol)

Description:

Ran (Ras-related nuclear protein, TC4) belongs to the superfamily of GTP binding proteins. It acts in cellular processes such as nuclear transport, mitotic spindle assembly, and post-mitotic nuclear assembly. The mutation T24N results in no or little binding to GTP and is known to act as a negative inhibitor of nuclear import. Protein preparation is 18% GDP-loaded, measured by HPLC.

Selected References:

Dasso (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.

Sekimoto *et al.* (1996) Interferon-gamma-dependent nuclear import of Stat1 is mediated by the GTPase activity of Ran/TC4. *J. Biol. Chem.* **271**:31017.