

RapGAP-341 (R91A)

human, recombinant, *E. coli*

Cat. No.	Amount
PR-222	50 µg

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

Avoid freeze / thaw cycles

Form

Liquid. Supplied in PBS and 5 mM DTE.

Molecular weight

39.5 kDa (amino acids 75 - 415)

Purity

> 95% by SDS-PAGE

Description

RapGAP, the GTPase activating protein of Rap, catalyzes the transition of Rap from GTP-bound to GDP-bound state by hydrolysis of GTP.

Rap proteins are GTPases of the Ras superfamily of G proteins opposing the Ras activation pathway. RapGap-341 is a catalytic active fragment comprising 341 amino acids (residues 75 - 415) of wild type Rap1GAP. RapGAP-341 (R91A) retained more than 50% activity in comparison to RapGAP-341.

Selected References:

Brinkmann et al. (2002) Rap-specific GTPase activating protein follows an alternative mechanism. *J. Biol. Chem.* **277**:12525.

Rubinfeld et al. (1992) Localization of the rap1GAP catalytic domain and sites of phosphorylation by mutational analysis. *Mol. Cell Biol.* **12**:4634.

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 GEFs, small GTPases, heterotrimeric G-proteins, and many more...

For detailed information please view the sections on

<http://www.jenabioscience.com>