

EPAC-1 Δ DEP

human, recombinant, *E. coli*

Cat. No.	Amount
PR-304	50 μ g

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

Avoid freeze / thaw cycles

Form

Liquid. Supplied in 50 mM Tris-HCl pH 7.6, 50 mM NaCl, 5 mM DTE and 5 % glycerol.

Molecular Weight

82.0 kDa (amino acids 149 - 883)

Purity

$\geq 95\%$ by SDS-PAGE

Description

EPAC (exchange protein directly activated by cAMP) is a Rap-specific guanine-nucleotide exchange factor (GEF). EPAC is activated by the binding of cAMP to a cyclic nucleotide monophosphate-binding domain.

N-terminal deletion of the DEP domain (domain that occur in Dishevelled, Egl-10, and Pleckstrin) of EPAC Δ DEP does not affect regulation of EPAC-activity but affects membrane localization.

Selected References:

Qiao *et al.* (2002) Cell cycle-dependent subcellular localization of exchange factor directly activated by cAMP. *J. Biol. Chem.* **277**:26581.

Krämer *et al.* (2001) Dynamic interaction of cAMP with the Rap guanine-nucleotide exchange factor EPAC1. *J. Mol. Biol.* **306**:1167.

De Rooij *et al.* (2000) Mechanism of regulation of the EPAC family of cAMP-dependent Rap GEFs. *J. Biol. Chem.* **275**:20829.

De Rooij *et al.* (1998) Epac is a Rap1 guanine-nucleotide-exchange factor directly activated by cAMP. *Nature* **396**:474.

Associated products available from Jena Bioscience

Non-hydrolyzable GTP-analogs (such as GTP γ 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.