

NORE1-RA^{GST}

Novel Ras effector 1, Ras Association Domain (AA 188-390)

human, recombinant, *E. coli*

Cat. No.	Amount
PR-307	200 µg

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

Avoid freeze / thaw cycles

Form

Liquid. Supplied in PBS supplemented with 1 mM EDTA.

Molecular weight

47 kDa with GST-tag.

Purity

>90% by SDS PAGE.

Description

Nore1 (Novel Ras effector 1) is a non-catalytic 413 amino acid polypeptide that contains several PXXP motifs, a central C1 zinc finger and a Ras-Rap association domain (RA) at its C-terminus.

Consistently, Nore1 interacts in a GTP-dependent fashion with a number of Ras-family GTPases, including Ras itself, M-Ras, TC21 and Rap2A (1,2).

The C-terminal portion of Nore1 exhibits 50 % identity with members of the RASSF Family of putative tumour suppressor proteins. Nore1 mediates the antiproliferative and/or pro-apoptotic effects of Ras in certain cell types, at least partly by activating the MST1 kinase (Mammalian Ste20-like kinase) (3).

Recent reports on the epigenetic silencing of Nore1 in a variety of human tumours have made a strong case for a role of Nore1 as a tumour suppressor (4,5).

Selected References:

- Ortiz-Vega, S. *et al.* (2002) *Oncogene*, **21**:1381.
Harjes, E. *et al.* (2006) *Structure*, **14**:881.
Praskova, M. *et al.* (2004) *Biochem J.*, **381**:453.
Tommasi, S.R. *et al.* (2002) *Oncogene*, **21**:2713.
Hesson, L. *et al.* (2003) *Oncogene*, **22**:947.