

## SRC1-RID<sup>GST</sup>

### Steroid Receptor Coactivator 1-Receptor Interacting Domain

human, recombinant, *E. coli*

Cat. No.	Amount
PR-831	10 µg

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

#### Avoid freeze / thaw cycles

#### Form

Liquid. Supplied in 20 mM Tris-HCl pH 8.0, 20% glycerol, 100 mM KCl, 0.2 mM EDTA and 1 mM DTT.

#### Activity

100 ng are sufficient for a protein-protein interaction assay.

#### Application

Use only for research and not for diagnostic purposes. Recombinant GST-SCR1 (627-786) can be used for protein-protein interaction assays.

#### Molecular Weight

58 kDa

#### Purity

> 95% by SDS-PAGE

#### Description

Steroid receptor coactivator 1 (SRC1) is a transcriptional coactivator that mediates the activating functions of many of the nuclear hormone receptors. It is also known as NCoA1 and is a member of the SRC/p160 coactivator family. SRC1 is a 160 kDa protein that contains several LXXLL motifs, which are involved in nuclear receptor interaction. The region 627-786 contains 3 LXXLL motifs that are involved in interaction with nuclear hormone receptors and has been previously used in assays detecting liganddependent receptor-cofactor interactions.

#### Selected References:

- Xu *et al.* (2003) Review of the *in vivo* functions of the p160 steroid receptor coactivator family. *Mol. Endocrinol.* **17**:1681.  
Bai *et al.* (2003) Isoform-selective interactions between estrogen receptors and steroid receptor coactivators promoted by estradiol and ErbB-2 signaling in living cells. *Mol. Endocrinol.* **17**:589.  
Cho *et al.* (2003) A simple ELISA for screening ligands of peroxisome proliferator-activated receptor-gamma. *J. Biochem. Mol. Biol.* **36**:207.  
Heery *et al.* (1997) A signature motif in transcriptional co-activators mediates binding to nuclear receptors. *Nature* **387**:733.  
Onate *et al.* (1995) Sequence and characterization of a coactivator for the steroid hormone receptor superfamily. *Science* **270**:1354.