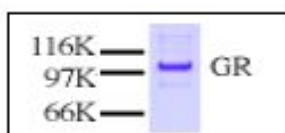


## GR

### Glucocorticoid Receptor

human, recombinant, Sf9 insect cells

Cat. No.	Amount
PR-750	5 $\mu$ g



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, 100 mM KCl, 0.2 mM EDTA, 1 mM DTT, 20% glycerol.

#### Activity

20 ng are sufficient for a gel-mobility shift assay and 100 ng are sufficient for a protein-protein interaction assay.

#### Application

GR has been applied in DNA and protein-protein interactions assays.

#### Purity

> 95% by SDS-PAGE.

#### Description

Recombinant His tagged GR was expressed in a baculovirus system and purified by affinity and FPLC chromatography.

Glucocorticoids are a vital class of steroid hormones that mediate profound and diverse physiological effects in vertebrates from fish to man. Although named for their role in glucose homeostasis, glucocorticoids are eminently important throughout physiology, with regulatory roles in development, metabolism, neurobiology, programmed cell death, and many other functions. In addition to these far-reaching physiological roles, corticosteroids are among the most widely prescribed class of drugs in the world. The physiological response and sensitivity to glucocorticoids varies among species, individuals, tissues, cell types, and even during the cell cycle. Additionally, several pathological conditions lead to, or are a result, of glucocorticoid resistance or hypersensitivity. The ligand-activated GR also interacts with a multitude of transcription factors such as c-jun, Nuclear Factor-B (NF-B), the TFIIID complex, STAT5, as well as a host of coactivators where they are known to act on the function of these signaling molecules. In addition, the GR interacts with numerous cytosolic proteins including chaperones, kinases, phosphatases, nuclear shuttling proteins, and the proteasome.

Recombinant GR was expressed in a baculovirus system and purified by an affinity column in combination with FPLC chromatography.

#### Selected References:

- Hsu *et al.* (1995) Selectivity of cell cycle regulation of glucocorticoid receptor function. *J. Biol. Chem.* **270**:3359.
- Lim-Tio *et al.* (1997) Determinants of specificity of transactivation by the mineralocorticoid or glucocorticoid receptor. *Endocrinology* **138**:2537.
- Kino and Chrousos (2001) Glucocorticoid and mineralocorticoid resistance/hypersensitivity syndromes. *Endocrinology* **169**:437.
- McKay and Cidlowski (1999) Molecular control of immune/inflammatory responses: interactions between nuclear factor-kappa B and steroid receptor-signaling pathways. *Endocr. Rev.* **20**:435.
- Ford *et al.* (1997) Involvement of the transcription factor IID protein complex in gene activation by the N-terminal transactivation domain of the glucocorticoid receptor *in vitro*. *Mol Endocrinol.* **11**:1467.
- Stocklin *et al.* (1996) Functional interactions between Stat5 and the glucocorticoid receptor. *Nature* **383**:726.
- Jenkins *et al.* (2001) Novel glucocorticoid receptor coactivator effector mechanisms. *Endocrinol. Metab.* **12**:122.
- Yudt and Cidlowski (2002) The glucocorticoid receptor: coding a diversity of proteins and responses through a single gene. *Mol. Endocrinol.* **16**:1719.