

## PPAR- $\beta$ -LBD (residues 165-441)

Peroxisome Proliferator Activated Receptor, beta isoform, Ligand Binding Domain  
human, recombinant, *E. coli*

| Cat. No. | Amount     |
|----------|------------|
| PR-846   | 10 $\mu$ 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, 20% glycerol, 100 mM KCl, 1 mM DTT and 0.2 mM EDTA.

#### Application

PPAR $\beta$  can be applied in DNA and protein-protein interactions assays.

#### Molecular Weight

30 kDa

#### Purity

> 95% by SDS-PAGE

### Description

Recombinant His tagged PPAR $\beta$ -LBD is isolated from an *E. coli* strain that carries the coding sequence of the human PPAR $\beta$  under the control of a T7 promoter.

There is evidence that a group of closely related nuclear receptors, called peroxisome proliferator-activated receptors (PPARs), may be involved in chronic diseases such as diabetes, obesity, atherosclerosis and cancer. The PPARs were first cloned as the nuclear receptors that mediate the effects of synthetic compounds called peroxisome proliferators on gene transcription. It soon became clear that eicosanoids and fatty acids can also regulate gene transcription through PPARs. They bind a specific element in the promoter region of target genes only as a heterodimer with the receptor for 9- cis retinoic acid, RXR (retinoid X receptor). Binding of the ligand of either receptor can activate the complex, but binding of both ligands simultaneously is more potent. Three PPAR isotypes have been identified:  $\alpha$ ,  $\beta$  (also called NUC1) and  $\gamma$ . PPAR $\alpha$  is expressed most in brown adipose tissue and liver, then kidney, heart and skeletal muscle.

PPAR $\gamma$  is mainly expressed in adipose tissue, and to a lesser extent in colon, the immune system and the retina. PPAR $\beta$  is found in many tissues but the highest expression is in the gut, kidney and heart. PPAR $\beta$  has received little attention, probably because of the lack of a connection with important clinical manifestations. However, recently PPAR $\beta$  has been linked to colon cancer, among other functions. PPAR regulates the expression of acyl-CoA synthetase 2 in the brain, linking PPAR $\beta$  to basic lipid metabolism. Moreover, it probably participates in embryo implantation and decidualization.

### Selected References:

- Kersten (2000) Roles of PPARs in health and disease. *Nature* **405**:421.
- Desvergne *et al.* (1999) Peroxisome proliferator-activated receptors: nuclear control of metabolism. *Endocr. Rev.* **20**:649.
- He *et al.* (1999) PPARdelta is an APC-regulated target of nonsteroidal anti-inflammatory drugs. *Cell* **99**:335.
- Basu-Modak *et al.* (1999) Peroxisome proliferator-activated receptor beta regulates acyl-CoA synthetase 2 in reaggregated rat brain cell cultures. *J. Biol. Chem.* **274**:35881.
- Lim *et al.* (1999) Cyclo-oxygenase-2-derived prostacyclin mediates embryo implantation in the mouse via PPARdelta. *Genes Dev.* **13**:1561.