

## PPAR- $\alpha$ -LBD (170-468)

**Peroxisome Proliferator-activated Receptor,  $\alpha$ -isoform**  
**human, recombinant, *E. coli***

Cat. No.	Amount
PR-816	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, 0.2 mM EDTA and 1 mM DTT.

#### Molecular Weight

33 kDa

#### Activity

20-100 ng are sufficient for a ligand binding assay and 100 ng are sufficient for a protein-protein interaction assay.

#### Application

PPAR- $\alpha$  has been applied in DNA and protein-protein interaction assays.

#### Purity

> 95% by SDS-PAGE

### Description

Recombinant His tagged PPAR $\alpha$ -LBD is isolated from an *E. coli* strain that carries the coding sequence of the human PPAR $\alpha$  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. The target genes of PPAR- $\alpha$  are a relatively homogenous group of genes that participate in aspects of lipid catabolism such as fatty acid uptake through membranes, fatty acid binding in cells, fatty acid oxidation (in microsomes, peroxisomes and mitochondria) and lipoprotein assembly and transport.

### Selected References:

- Kersten *et al.* (2001) Mechanisms of nutritional and hormonal regulation of lipogenesis. *EMBO Rep.* **21**:282.  
Kersten *et al.* (2000) Roles of PPARs in health and disease. *Nature* **405**:421.