

## Recombinant PXR protein

---

**Catalog No:** 31144

**Expressed In:** Baculovirus

**Quantity:** 5 µg

**Concentration:** 0.65 µg/µl

**Source:** Human

**Buffer Contents:** 5 µg of Recombinant PXR protein in Dilution Buffer AM1 (20 mM Tris-Cl (pH 7.5), 20% glycerol, 100 mM KCl, 1 mM DTT and 0.2 mM EDTA).

**Protein Details:** Recombinant PXR (accession number NM\_003889) is expressed in a baculovirus system; carries the full-length coding sequence and is purified by an affinity column in combination with FPLC chromatography.

**Application Notes:** PXR can be used for Western blot (50 kDa), gel mobility shift assay, for protein-protein and small molecule-protein interaction assays. 1 ng is sufficient for a gel mobility shift assay in a 20 µl reaction; 100 ng is sufficient for a protein-protein interaction assay detected by immuno-blot system.

NOTE: The presence of Poly [d(I-C)] in buffers may affect protein functionality and should be avoided.

### References:

This product was used in the following publications:

Wang, C., *et. al.* (2018). "Poly(ADP-ribosyl)ated PXR is a critical regulator of acetaminophen-induced hepatotoxicity." *Cell Death Dis.* 9(8):819. PMID: 30050067. (Western Blot).

**Storage and Guarantee:** Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.