

# TREVIGEN<sup>®</sup> Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

## Recombinant Peripheral benzodiazepine receptor (PBR) protein

**Catalog#:** 6360-025-01

**Size:** 25  $\mu$ l

**Description:** Recombinant full length mouse peripheral benzodiazepine receptor (PBR) purified from *E. coli*.

**Applications:** Control for Western Blot analysis using the anti-PBR antibody.

**Physical State:** This purified protein is provided at 0.2 mg/ml in 20 mM Tris 7.8, 1 mM EDTA, 0.1% Sodium Lauryl Sarcosine.

**Expected Size:** The recombinant protein is a fusion protein with a molecular weight of 20.4 kDa. Please note that multimers of the proteins are commonly seen on gels and Western Blotting Detection. The native protein is a 18 kDa protein.

**Storage:** Store at -20°C in working aliquots in a manual defrost freezer to avoid repeated freeze-thaws.

### To use as a Western blot control:

Use 1  $\mu$ g of the recombinant PBR protein per lane (5 $\mu$ l). Dilute the protein with 1X loading buffer of choice, supplemented with 2 mM DTT. Heat denature the control protein for 5 minutes before loading.

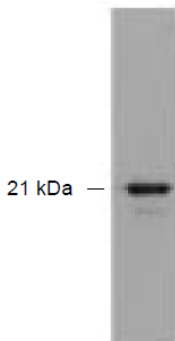


Fig.1 Western Blot analysis of PBR where 5  $\mu$ l of the recombinant PBR protein was loaded on a 12% SDS-PAGE gel, transferred to a PVDF membrane and detected with a 1:1000 dilution of the anti-PBR antibody (cat# 6362-PC-100). Detection was performed with a peroxidase-conjugated secondary antibody to rabbit followed by chemiluminescence.

## TREVIGEN<sup>®</sup>

8405 Helgerman Court, Gaithersburg, MD 20877 USA

Voice: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800

Fax: 301-560-4973 • e-mail: [info@trevigen.com](mailto:info@trevigen.com) • [www.trevigen.com](http://www.trevigen.com)

**Recombinant  
PBR Protein**  
Catalog #: 6360-025-01  
Storage: -20 °C

**TREVIGEN®**  
1-800-873-8443