

Data Sheet

FXR, GST-Tag

Human, Recombinant, N-terminal GST-tag
Catalog #: 100410
Lot #: 190717-1 **Conc.:** 1.27 mg/ml

Description: Human FXR also known as NR1H4, Bile Acid Receptor, and HRR1, GenBank Accession No. NM_005123, a.a. 193-472(end), with N-terminal GST-tag, expressed in a baculovirus infected Sf9 cell expression system. MW=59 kDa.

Formulated in: 50 mM Tris, pH 8, 100mM KCl, 5 mM DTT, 0.5 mM EDTA, 20% Glycerol, 1mM Glutathione.

Stability: At least 6 months at -80°C. Avoid freeze/thaw cycles.

References:

1. Cai, S.-Y., *et al.*, *Comp. and Evol. Phys.* 2007 Sep 1;**293(3)**:R1400-R1409.
2. Attinkara, R. *et al.*, *BMC.* 2012 Aug 28;**5**:461.

Quality Assurance

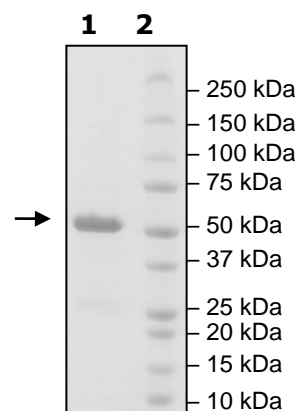
4-20% SDS-PAGE Coomassie staining

Lane 1:
FXR

Lane 2:
Protein Marker

Purity: ≥90%

MW: 59 kDa



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com