

Data Sheet

LDLR, FLAG-Tag

Human, Recombinant, C-terminal FLAG-tag

Catalog #: 71205

Lot #: 170223-2 **Conc.:** 1.10 mg/ml

Formulated in: 8 mM Phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol, 100µg/mL FLAG peptide.

Stability: At least 6 months at -80°C. Avoid freeze thaw cycles. Storing diluted protein is not recommended, if necessary, use a carrier protein (BSA 0.1 – 0.5%).

References:

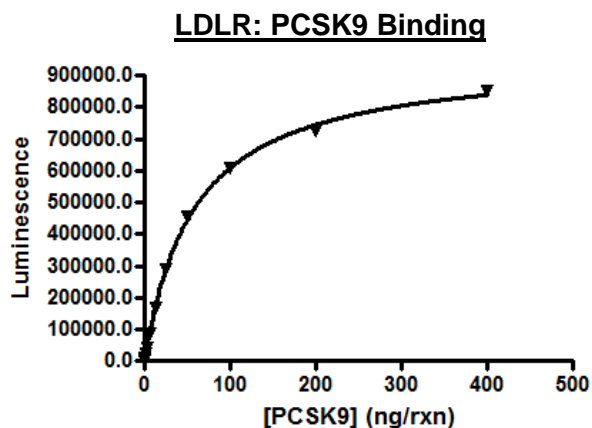
1. Holla, L., *et al.*, *BMC Cell Biol.* 2007 Mar 1;**8**:9.
2. Qian, YW., *et al.*, *J Lipid Res.* 2007 Jul;**48**(7):1488-98.
3. Fasano, T., *et al.*, *Athersclerosis.* 2009 Mar;**203**(1):166-71.

Description: Human low density lipoprotein receptor (LDLR), also known as FH, FHC, and LDLCQ2, GenBank Accession No. NM_000527, a.a. 22-788, with C-terminal Flag-tag, MW=86 kDa, expressed in a HEK293 cell expression system. LDLR is heavily glycosylated, resulting in higher molecular weight. The two bands shown correspond to differing states of glycosylation.

Assay Conditions: 100 ng/well of LDLR in 50 µl (overnight at 4°C). Initiated binding reaction with addition of various concentrations of PCSK9-biotin (50 ng/well is the established kit condition). Incubated for 2 hours at room temperature. Detected binding with Strep-HRP generated luminescence.

Application: Useful for studying protein binding and screening small molecules.

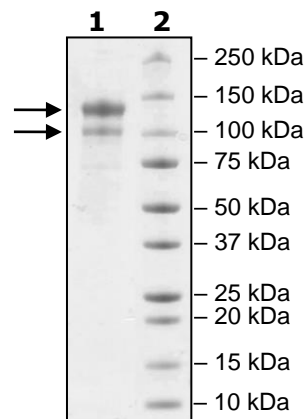
Quality Assurance



**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
2 µg LDLR
Lane 2:
Protein Marker

MW: 86 kDa +
Glycans
Purity: ≥90%



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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com