

Data Sheet

LDLR, FLAG-Avi-Tag Biotin-Labeled

Human, Recombinant, C-terminal Flag-Avi-tags
 Biotin-Labeled

Catalog #: 71206

Lot #: 190806-B **Conc.:** 0.71 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
 110 mM NaCl, 2.2 mM KCl, and 20%
 glycerol.

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

References:

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

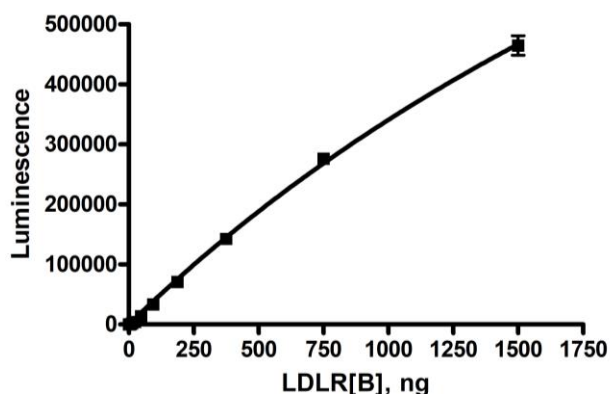
Description: Human low density lipoprotein
 receptor (LDLR), also known as FH, FHC, and
 LDLCQ2, GenBank Accession No. NM_000527,
 a.a. 22-788(end), with C-terminal FLAG and Avi
 tags, MW=88 kDa*, expressed in a HEK293 cell
 expression system and enzymatically
 biotinylated using Avi-tag™ technology. *Note:
 LDLR is heavily glycosylated, resulting in higher
 molecular weight. The two bands shown
 correspond to differing states of glycosylation.

Assay Conditions: Coat a plate with 100 ng of
 PCSK9. Add a 25 μL reaction mix (50 mM
 HEPES pH 7.4, 50 mM NaCl, 200 μM CaCl₂,
 0.05% Tween, 0.1% BSA, and up to 200 ng
 LDLR-biotin) to the coated wells and incubate
 for 2 hr. Add Streptavidin-HRP labeled
 antibody and incubate 1 hr. Finally, add HRP
 chemiluminescent substrates and read
 luminescence.

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

Quality Assurance

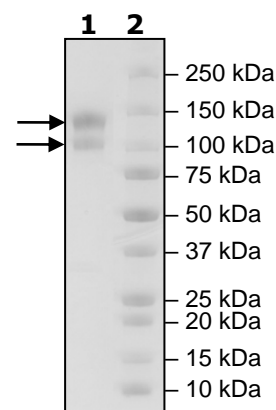
PCSK9: LDLR [Biotin] Interaction



4-20% SDS-PAGE Coomassie staining

Lane 1:
 4 μg LDLR
Lane 2:
 Protein Marker

MW: 88 kDa+glycans
Purity: $\geq 90\%$



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