

Data Sheet

LDLR, FLAG-Avi-Tag (Mouse)

Mouse, Recombinant, C-Terminal
FLAG-tag, C-terminal Avi-Tag

Catalog #: 71222

Lot #: 130110-D **Conc.:** 0.1 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,
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.

Description: Mouse low density lipoprotein
receptor (LDLR), also known as FH, FHC,
and LDLCQ2, GenBank Accession No.
NM_010700, a.a. 22-790, with C-terminal
Flag-tag, MW=87 kDa, expressed in a
HEK293 cell expression system. *Note:
LDLR is heavily glycosylated, resulting in
higher molecular weight.

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

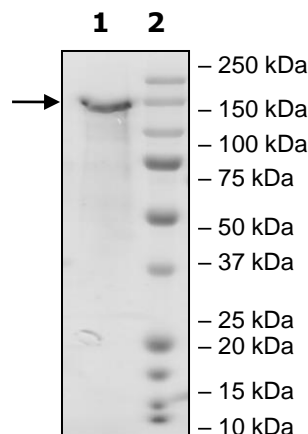
Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
LDLR

Lane 2:
Protein Marker

MW: 87 kDa+glycans
Purity: $\geq 88\%$



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