

Data Sheet

FNDC5, Avi-His-Tag, Biotin-Labeled

Human, recombinant, C-terminal Avi-His-Tag
Biotin labeled

Catalog #: 100426

Lot #: 190711-1

Conc.: 0.79 mg/ml

Formulated in: 8 mM phosphate, pH 7.4,
110 mM NaCl, 2.2 mM KCl, and 20%
glycerol.

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

References:

1. Wrann, C., *et al.* 2013. *Cell Metabolism*.
18(5): e60563
2. Roca-Rivada, A., *et al.* 2013. *J. PLoS*
ONE. **8(4):** 29-32.

Description: Human recombinant
Fibronectin type III domain-containing
protein 5 (FNDC5), also known as Irisin,
Fibronectin type III repeat-containing protein
2, and FRCP2, GenBank Accession No.
NM_153756, a.a 32-143, with a C-terminal
Avi-His-Tag, expressed in a HEK293 cell
expression system and enzymatically biotin-
labeled using Avi-tag™ technology.
Biotinylation is confirmed to be $\geq 90\%$.
MW=16 kDa. This protein runs at a higher
MW by SDS-PAGE due to glycosylation.

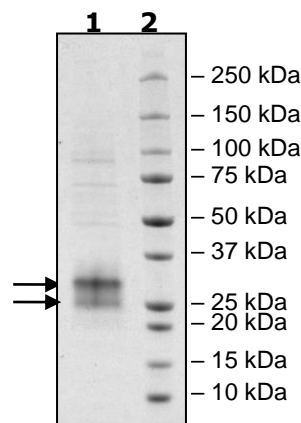
Quality Assurance

4-20% SDS-PAGE
Coomassie staining

Lane 1:
5 μg FNDC5

Lane 2:
Protein Marker

MW: 16 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 **1.858.202.1401** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com