

Data Sheet

FNDC5, Avi-His-Tag

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

Catalog #: 100425

Lot #: 190709-1

Conc.: 0.19 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. MW=16 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Quality Assurance

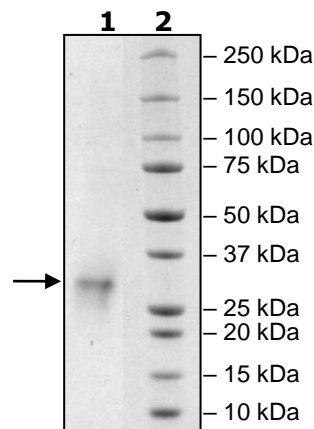
4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg FNDC5

Lane 2:
Protein Marker

MW: 16 kDa

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