

PRODUCT INFORMATION



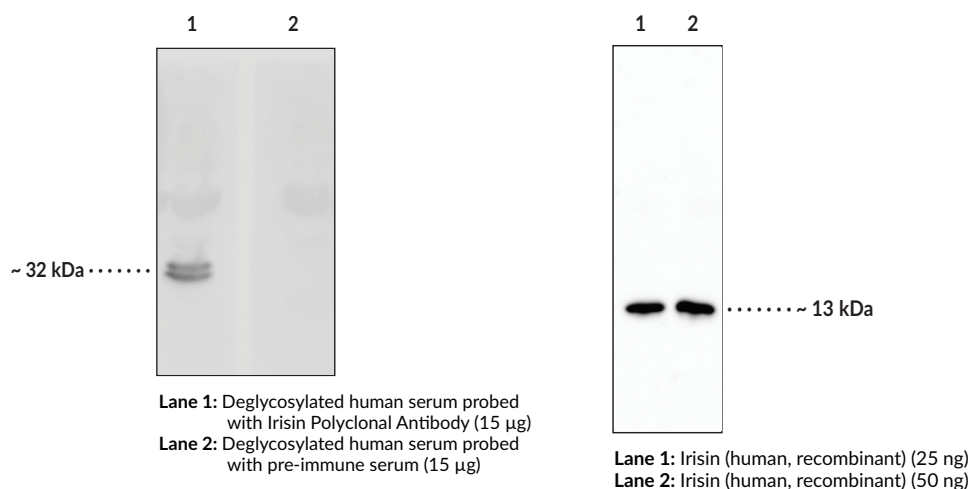
Irisin Polyclonal Antibody

Item No. 14625

Overview and Properties

Contents: This vial contains protein A-purified polyclonal antibody.
Synonyms: Fibronectin type III domain-containing protein 5, FNDC5, FRCP2, PeP
Immunogen: Synthetic peptide from within human irisin
Species Reactivity: (+) Human; other species not tested
Uniprot No.: Q8NAU1
Form: Lyophilized
Storage: -20°C (as supplied)
Stability: ≥2 years
Storage Buffer: TBS, pH 7.4 when reconstituted in 500 µl deionized water
Host: Rabbit
Applications: ELISA and Western blot (WB); the recommended starting dilution for ELISA is 1:500-1:2,000 and WB is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/04/2021

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Physical exertion and exercise are the primary defenses against obesity, insulin resistance, and diabetes. The transcriptional co-activator peroxisome proliferator-activated receptor C coactivator 1 α (PGC-1 α), regulates mitochondrial biogenesis and function. Expression of the membrane protein fibronectin type III domain-containing protein 5 (FNDC5) is stimulated in muscle by increased expression of the PGC1- α protein in response to exercise. FNDC5 is proteolytically cleaved and secreted as the hormone peptide irisin (named after the Greek goddess messenger Iris). Irisin has been shown to convert white adipose tissue (WAT) to brown adipose tissue (BAT).¹ An increase in BAT results in anti-obesity and anti-diabetic metabolic states in mice. Exogenous irisin levels in mice leads to higher levels of the uncoupling protein UCP1, which results in an increase in thermogenesis and heat expenditure.² The physiological responses induced by the effects of irisin have the potential to increase weight loss and reduce insulin resistance and metabolic disorder. Cayman's irisin polyclonal antibody detects a 13 kDa band from bacterially expressed irisin protein (Item No. 11451), and a 32 kDa band from deglycosylated human serum.

References

1. Boström, P., Wu, J., Jedrychowski, M.P., *et al.* A PGC1- α -dependent myokine that drives brown-fat-like development of white fat and thermogenesis. *Nature* **481**, 463-468 (2012).
2. Wu, J., Boström, P., Sparks, L.M., *et al.* Beige adipocytes are a distinct type of thermogenic fat cell in mouse and human. *Cell* **150**, 366-376 (2012).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM