

Irisin (human) polyclonal antibody

Irisin is a hormone that is involved in converting white adipose tissue into brown adipose tissue after physical exercise. It is formed after the membrane bound protein fibronectin type III domain containing protein 5 (FNDC5) is cleaved resulting in a secreted protein. FNDC5 is found in muscle tissue and is activated during physical exertion via the protein PGC1a. Soluble Irisin has also been shown to up regulate UCP-1 gene expression in mice. UCP-1 is a mitochondrial transporter protein and a marker of brown adipose tissue conversion. This suggests that Irisin plays an important role in energy regulation and furthermore may lead to improvements in glucose regulation, insulin sensitivity, hypertension, and obesity.

This antibody is covered by our [Worry-Free Guarantee](#).

Ordering Information

[Order Online »](#)

ENZ-ABS256-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Avoid freeze/thaw cycles.

Short Term Storage -20°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	FNDC5, Fibronectin type III domain-containing protein 5, FRCP2, Fibronectin type III repeat-containing protein 3
Application	ELISA, IP, WB
Formulation	Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.
Host	Rabbit
Immunogen	Recombinant human Irisin (active).
Positive Control	ADI-908-307
Purity Detail	Protein A affinity purified and immunodepleted.
Source	Purified from rabbit serum.
Species Reactivity	Human
UniProt ID	Q8NAU1
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee



ENZO LIFE SCIENCES,
Last modified **May 29, 2024**
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-be@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com