

IGF-I (human), (recombinant)

Recombinant human IGF-I that promotes cell proliferation and inhibits apoptosis.

Insulin-like growth factor-1 (IGF-1) was originally known as Somatomedin C but was later changed to IGF-1 due to its structural similarity to proinsulin. IGF-1 has insulinlike effects, but it appears to be involved in many other functions and has been linked to aging, nutrition, physical activity and cancer. While each of these is very interesting on its own, together they may shed light on the role that lifestyle plays in influencing cancer risk. High levels of circulating IGF-1 and low levels of IGF binding protein-3 (IGFBP-3) are associated with increased risk of several cancers. IGF-1 can stimulate cell proliferation and inhibit apoptosis in many different kinds of cells. IGF-1 circulates in blood bound with IGFBPs. Six IGFBPs have been identified but the majority of circulating IGF-1 is complexed with IGFBP-3. The IGF-1 and IGFBP-3 exist in a complex with a third protein, named the acid-labile subunit.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ADI-908-059-0100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Lyophilized IGF-I although stable at room temperature for 3 weeks, should be stored desiccated below -18°C. Upon reconstitution IGF-I should be stored at +4°C between 2-7 days and for future use below -18°C. For long term storage it is recommended to add a carrier protein (0.1% HSA or BSA).
Handling	Avoid freeze-thaw cycles.
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Insulin-like growth factor I
Application Notes	Western blot control.
Biological Activity	The activity as determined by the dose-dependent proliferation of mouse FDC-P1 is less than 1.0 ng/ml, corresponding to a specific activity of 1×10^6 units/mg.
Formulation	Lyophilized from a sterile filtered concentrated solution in PBS, pH 7.0.
MW	~8kDa
Purity	≥95% (SDS and HPLC)
Purity Detail	Purified by SDS and HPLC.
Reconstitution	Reconstitute the lyophilized IGF-1 in sterile 18MΩ-cm H ₂ O not less than 100µg/ml, which can then be further diluted to other aqueous solutions.
Source	Produced in <i>E. coli</i> .

UniProt ID



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

P05019

France
Phone: +33 472 440 655
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

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

Last modified: October 9, 2025

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.